

# PARP-1 monoclonal antibody (C-2-10) (affinity purified)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 36

[View Online »](#)

## Ordering Information

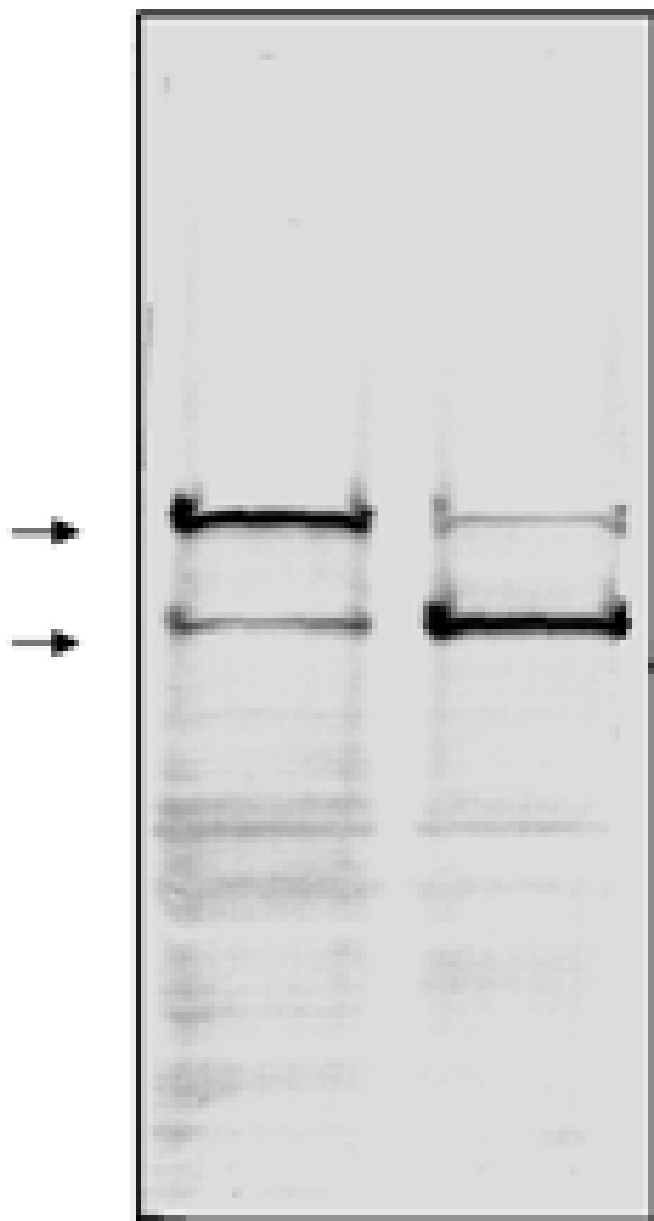
[Order Online »](#)

BML-SA249-0050	50µg
----------------	------

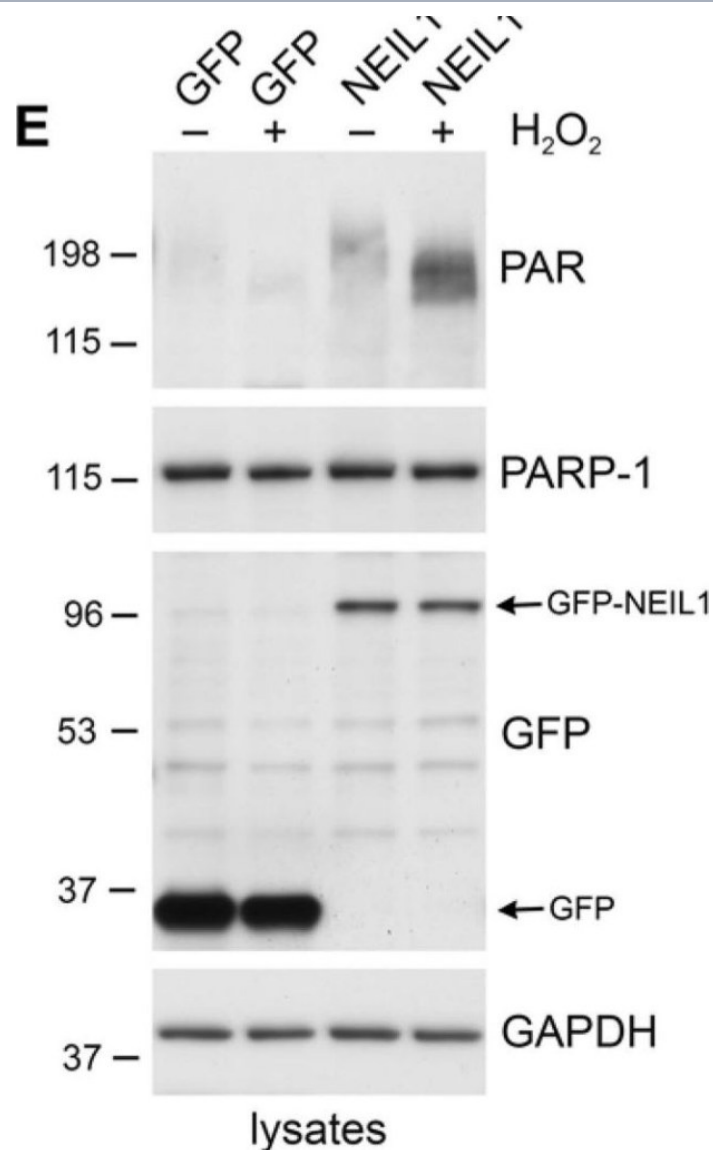
Manuals, SDS & CofA

[View Online »](#)

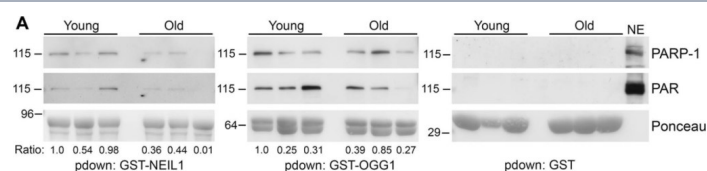




Western blot analysis of whole cell extracts from (1) human HL60 leukemia cells (Prod. No. BML-SW101) and (2) HL60 cells induced to undergo apoptosis using etoposide (Prod. No. BML-SW102). The nitrocellulose membrane was probed with purified clone C-2-10 mouse monoclonal anti-PARP (Prod. No. BML-SA249) at 1  $\mu$ g/ml. Secondary antibody was GAM-AP (1/2000) and the nitrocellulose membrane was developed with BCIP/NBT. Arrows correspond to 116kDa and 85kDa.

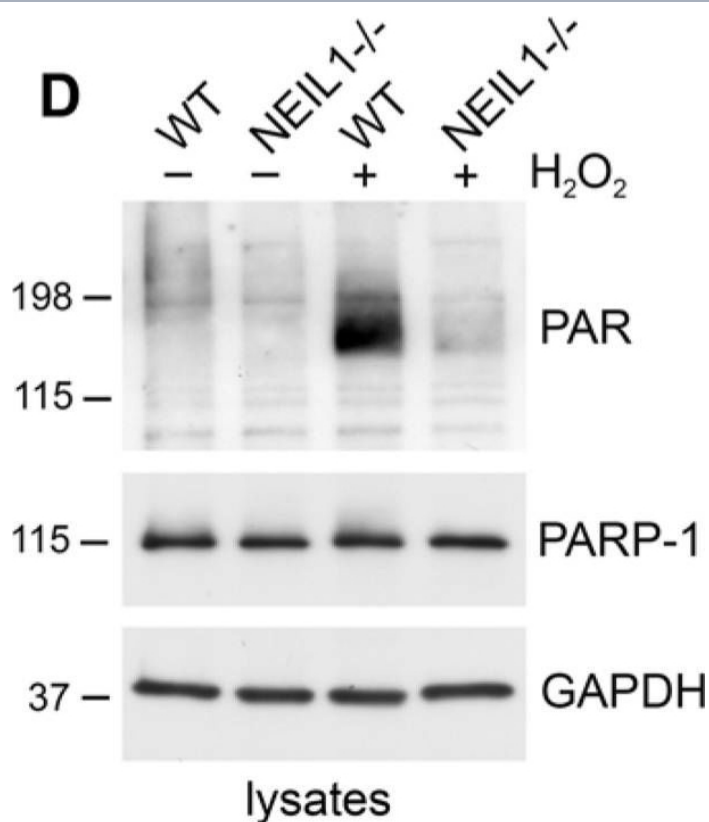


NEIL1 stimulates the poly(ADP-ribosyl)ation activity of PARP-1(A) PARP-1 activity was measured by determining the amount of PAR deposited on immobilized histones in an ELISA assay. Addition of NEIL1 (33 ng) increased the amount of PAR synthesis by PARP-1 (2 ng). The histogram represents the mean + SEM from three independent experiments. \* $p < 0.05$  compared to control by Student's t-test. (B,C) GST, GST-NEIL1, GST-NEIL1 1–288 aa or GST-NEIL1 289–390 aa (5  $\mu$ g) were incubated with (+) or without (–) 7.5 ng PARP-1. The immunoblot was stained with Ponceau S and probed with anti-PAR antibodies. (D) WT or NEIL1 $^{-/-}$  MEFs were mock treated (–) or treated (+) with 500  $\mu$ M H<sub>2</sub>O<sub>2</sub>. Lysates were probed with anti-PAR antibodies and reprobed with anti-PARP-1 and anti-GAPDH antibodies as a protein loading control. (E) NEIL1 $^{-/-}$  MEFs transfected with pEGFP-NEIL1 or pEGFP control were mock treated (–) or treated (+) with 500  $\mu$ M H<sub>2</sub>O<sub>2</sub> and lysates were probed as above. Anti-GFP antibodies were used to examine levels of the transfected proteins. For (B-E), a representative experiment from at least 3 independent experiments is shown.



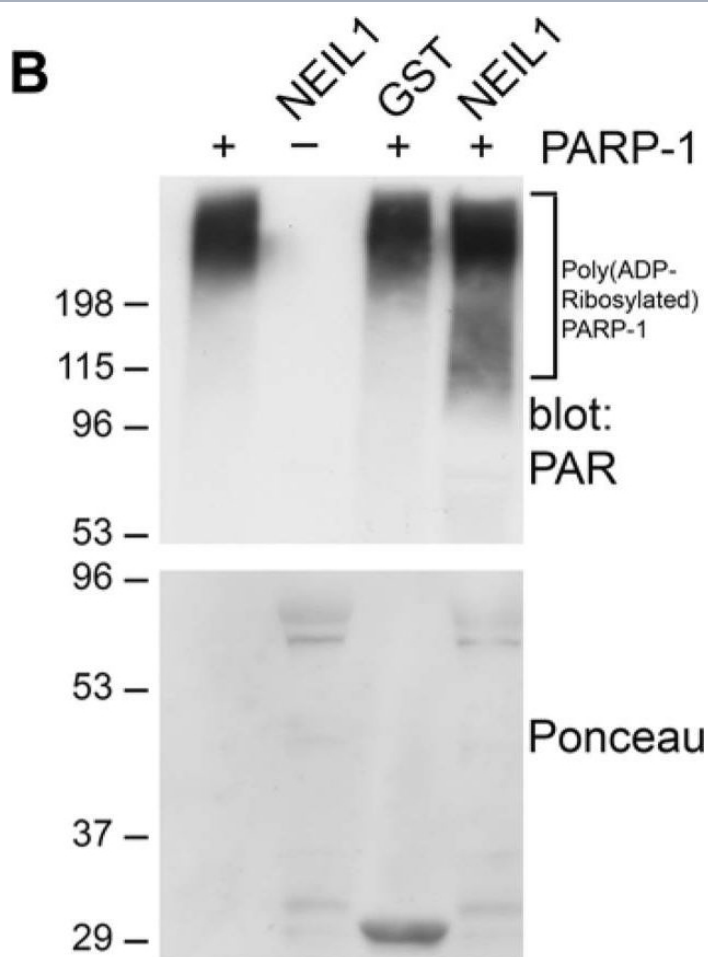
Decreased binding of PARP-1 to GST-NEIL1 with mouse age(A) Liver nuclear extracts were made from 3 different young (7 month) and 3 different old (17 month) mice and used for GST-precipitations with the indicated fusion proteins. Precipitations were probed with anti-PARP-1 and anti-PAR antibodies and stained with Ponceau S to visualize the fusion proteins. The relative amount of PARP-1 binding and PAR levels in precipitations were quantified from immunoblots and normalized to the amount of GST-fusion protein. The numbers below the blots represent the relative level of PARP-1 binding and the histograms in (B) represent the average PARP-1 binding and PAR levels in precipitations from the 3 young and 3 old mice. \*P<0.05 comparing young and old mice using Student's t-test. Similar results were obtained in another independent experiment. (C) Liver nuclear extracts from young and old mice were probed by immunoblotting with anti-PARP-1 and anti-NEIL1 antibodies and reprobed with anti-TBP antibodies as a protein loading control. The numbers below the blots represent the level of PARP-1 or NEIL1 normalized to TBP. The histogram in (D) represents the average levels of NEIL1 and PARP-1 + SEM from the 3 young and 3 old mice.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Coordination of DNA repair by NEIL1 and PARP-1: a possible link to aging. *Aging (Albany NY)* (2012)



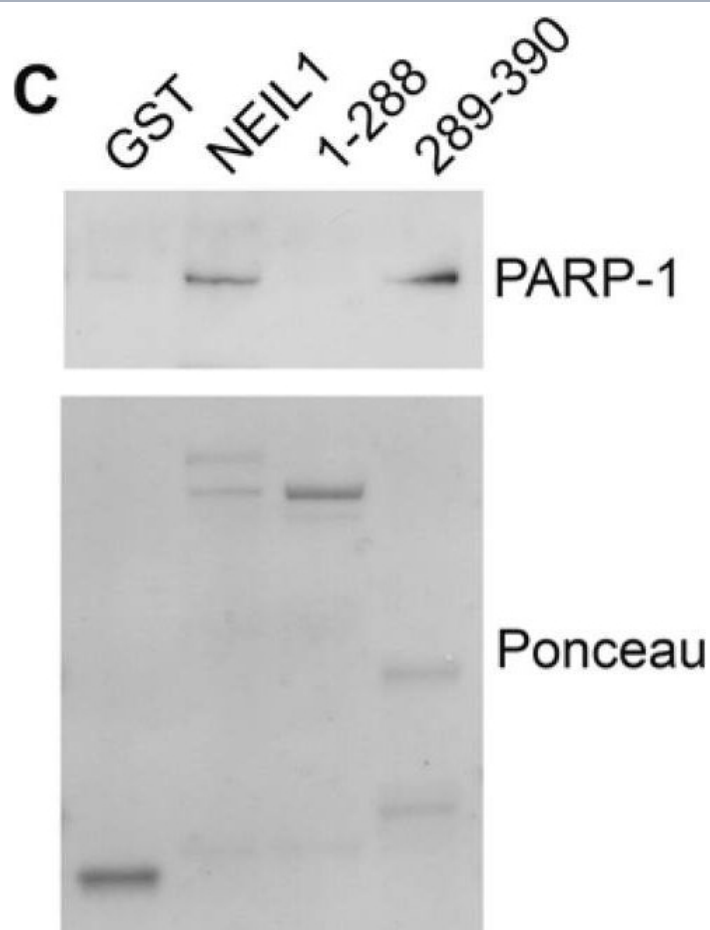
NEIL1 stimulates the poly(ADP-ribosyl)ation activity of PARP-1(A) PARP-1 activity was measured by determining the amount of PAR deposited on immobilized histones in an ELISA assay. Addition of NEIL1 (33 ng) increased the amount of PAR synthesis by PARP-1 (2 ng). The histogram represents the mean + SEM from three independent experiments. \*p < 0.05 compared to control by Student's t-test. (B,C) GST, GST-NEIL1, GST-NEIL1 1–288 aa or GST-NEIL1 289–390 aa (5 µg) were incubated with (+) or without (–) 7.5 ng PARP-1. The immunoblot was stained with Ponceau S and probed with anti-PAR antibodies. (D) WT or NEIL1–/– MEFs were mock treated (–) or treated (+) with 500 µM H<sub>2</sub>O<sub>2</sub>. Lysates were probed with anti-PAR antibodies and reprobed with anti-PARP-1 and anti-GAPDH antibodies as a protein loading control. (E) NEIL1–/– MEFs transfected with pEGFP-NEIL1 or pEGFP control were mock treated (–) or treated (+) with 500 µM H<sub>2</sub>O<sub>2</sub> and lysates were probed as above. Anti-GFP antibodies were used to examine levels of the transfected proteins. For (B–E), a representative experiment from at least 3 independent experiments is shown.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Coordination of DNA repair by NEIL1 and PARP-1: a possible link to aging. *Aging (Albany NY)* (2012)



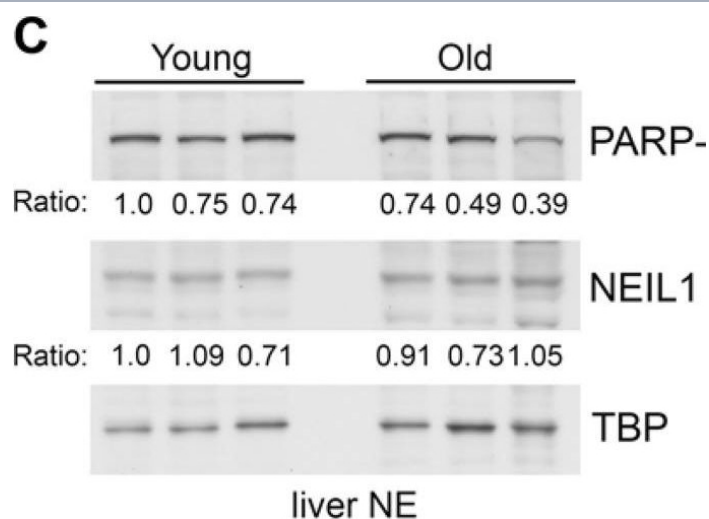
NEIL1 stimulates the poly(ADP-ribosylation) activity of PARP-1(A) PARP-1 activity was measured by determining the amount of PAR deposited on immobilized histones in an ELISA assay. Addition of NEIL1 (33 ng) increased the amount of PAR synthesis by PARP-1 (2 ng). The histogram represents the mean + SEM from three independent experiments. \* $p < 0.05$  compared to control by Student's t-test. (B,C) GST, GST-NEIL1, GST-NEIL1 1–288 aa or GST-NEIL1 289–390 aa (5  $\mu$ g) were incubated with (+) or without (–) 7.5 ng PARP-1. The immunoblot was stained with Ponceau S and probed with anti-PAR antibodies. (D) WT or NEIL1 $^{-/-}$  MEFs were mock treated (–) or treated (+) with 500  $\mu$ M H<sub>2</sub>O<sub>2</sub>. Lysates were probed with anti-PAR antibodies and reprobed with anti-PARP-1 and anti-GAPDH antibodies as a protein loading control. (E) NEIL1 $^{-/-}$  MEFs transfected with pEGFP-NEIL1 or pEGFP control were mock treated (–) or treated (+) with 500  $\mu$ M H<sub>2</sub>O<sub>2</sub> and lysates were probed as above. Anti-GFP antibodies were used to examine levels of the transfected proteins. For (B–E), a representative experiment from at least 3 independent experiments is shown.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Coordination of DNA repair by NEIL1 and PARP-1: a possible link to aging. *Aging (Albany NY)* (2012)



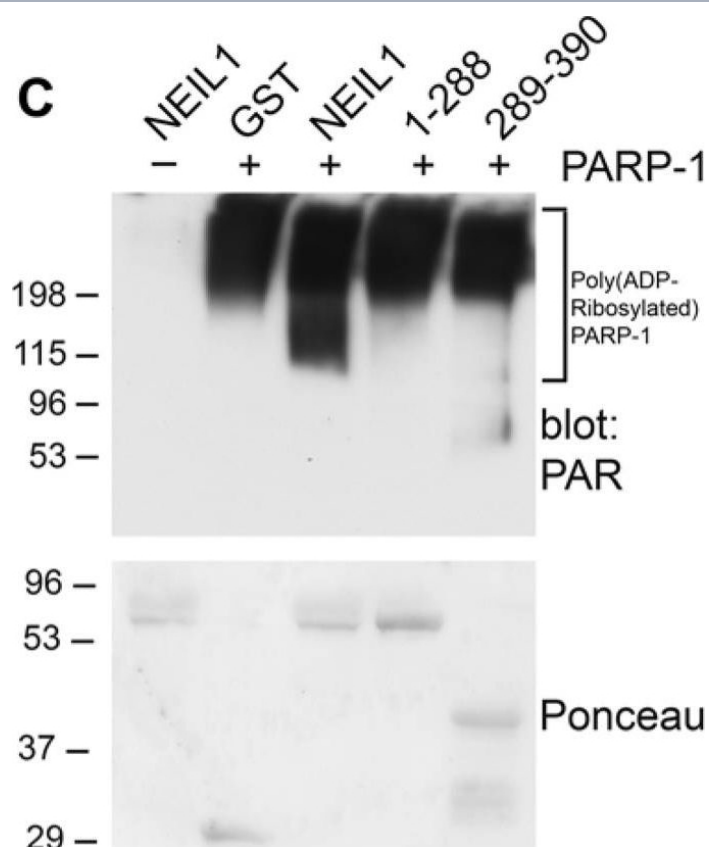
Mapping the interacting regions of PARP-1 and NEIL1(A) Schematic representation of GST-NEIL1 fusion proteins. (B) Purified GST-NEIL1 fusion proteins were analyzed by SDS-PAGE followed by staining with Coomassie Colloidal Blue. (C) PARP-1 binds to the C-terminal 289–390 aa of NEIL1. PARP-1 (250 ng) was incubated with either GST, GST tagged full-length NEIL1 or indicated NEIL1 fragments (1  $\mu$ g) in an in-vitro binding assay. GST precipitations were probed with anti-PARP-1 antibodies and stained with Ponceau S. ~5% of purified PARP-1 bound to both GST-NEIL and NEIL 289–390aa. (D) NEIL1 binds to the BRCT domain of PARP-1. The DNA binding (41 kDa), BRCT (15 kDa) and catalytic (39 kDa) domains of PARP-1 (1  $\mu$ g) were incubated with 1  $\mu$ g GST-NEIL1 (69 kDa) either in the presence (+) or absence (–) of ethidium bromide. GST-NEIL1 precipitations were immunoblotted with anti-His (for the DNA binding and catalytic domains) and anti-BRCT (for BRCT domain) antibodies and then stained with Coomassie to reveal the amount of GST-NEIL1 in the precipitations. The different PARP-1 domains and GST-NEIL1 are indicated by arrows. ~4% of purified BRCT domain bound to GST-NEIL1. (E) A schematic of the different PARP-1 proteins is shown.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Coordination of DNA repair by NEIL1 and PARP-1: a possible link to aging. *Aging (Albany NY)* (2012)



Decreased binding of PARP-1 to GST-NEIL1 with mouse age(A) Liver nuclear extracts were made from 3 different young (7 month) and 3 different old (17 month) mice and used for GST-precipitations with the indicated fusion proteins. Precipitations were probed with anti-PARP-1 and anti-PAR antibodies and stained with Ponceau S to visualize the fusion proteins. The relative amount of PARP-1 binding and PAR levels in precipitations were quantified from immunoblots and normalized to the amount of GST-fusion protein. The numbers below the blots represent the relative level of PARP-1 binding and the histograms in (B) represent the average PARP-1 binding and PAR levels in precipitations from the 3 young and 3 old mice. \* $P < 0.05$  comparing young and old mice using Student's t-test. Similar results were obtained in another independent experiment. (C) Liver nuclear extracts from young and old mice were probed by immunoblotting with anti-PARP-1 and anti-NEIL1 antibodies and reprobed with anti-TBP antibodies as a protein loading control. The numbers below the blots represent the level of PARP-1 or NEIL1 normalized to TBP. The histogram in (D) represents the average levels of NEIL1 and PARP-1 + SEM from the 3 young and 3 old mice.

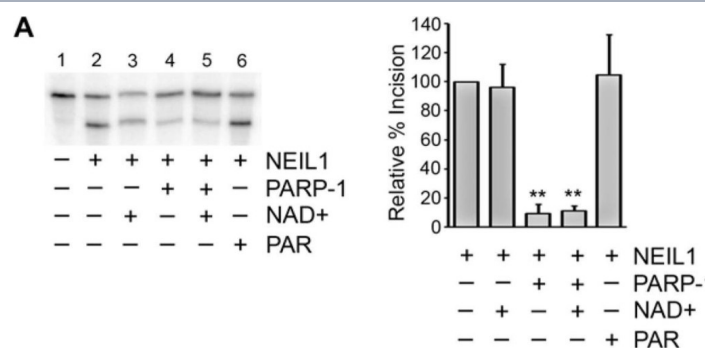
Image collected and cropped by CiteAb under a CC-BY license from the following publication: Coordination of DNA repair by NEIL1 and PARP-1: a possible link to aging. *Aging (Albany NY)* (2012)



NEIL1 stimulates the poly(ADP-ribosyl)ation activity of PARP-1(A) PARP-1 activity was measured by determining the amount of PAR deposited on immobilized histones in an ELISA assay. Addition of NEIL1 (33 ng) increased the amount of PAR synthesis by PARP-1 (2 ng). The histogram represents the mean + SEM from three independent experiments. \* $p < 0.05$  compared to control by Student's t-test. (B,C) GST, GST-NEIL1, GST-NEIL1 1–288 aa or GST-NEIL1 289–390 aa (5  $\mu$ g) were incubated with (+) or without (–) 7.5 ng PARP-1. The immunoblot was stained with Ponceau S and probed with anti-PAR antibodies. (D) WT or NEIL1 $^{-/-}$  MEFs were mock treated (–) or treated (+) with 500  $\mu$ M H<sub>2</sub>O<sub>2</sub>. Lysates were probed with anti-PAR antibodies and reprobed with anti-PARP-1 and anti-GAPDH antibodies as a protein loading control. (E) NEIL1 $^{-/-}$  MEFs transfected with pEGFP-NEIL1 or pEGFP control were mock treated (–) or treated (+) with 500  $\mu$ M H<sub>2</sub>O<sub>2</sub> and lysates were probed as above. Anti-GFP antibodies were used to examine levels of the transfected proteins. For (B–E), a representative experiment from at least 3 independent experiments is shown.

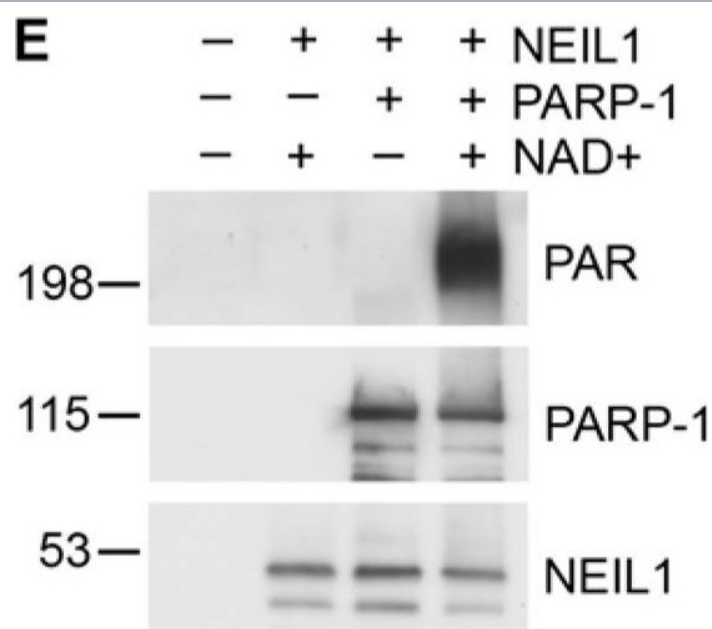
Image collected and cropped by CiteAb under a CC-BY license from the following publication: Coordination of DNA repair by NEIL1 and PARP-1: a possible link to aging. *Aging (Albany NY)* (2012)





PARP-1 inhibits NEIL1 incision activity(A) NEIL1 (38 nM; 33 ng) was incubated with buffer, NAD<sup>+</sup> with or without PARP-1 (180 nM; 400 ng), or with PAR and reacted with a 5'- 32P-labeled oligonucleotide duplex containing a 5OHU lesion for 15 min at 37°C. The cleavage products were analyzed on a 20% denaturing gel containing 7 M urea. Percent incision was calculated by normalizing the amount of cleaved substrate (bottom band) to the amount of uncleaved product (top band). Data was then normalized to the amount of incision activity of NEIL1 alone (100%). (B) Incision assays were performed as in (A) in the presence of the indicated concentrations of PARP-1 with (activated PARP-1) or without NAD<sup>+</sup> (PARP-1) and quantified. (C) NEIL1 (33 ng) incision assays were performed in the presence of 400 ng of the indicated PARP-1 domains or full-length PARP-1 in the presence of NAD<sup>+</sup> (10 μM). (D) PARP-1 inhibits NEIL1 activity at both 5 nM and 20 nM concentrations of DNA substrate. The histograms in (A) and (B) represent the mean + SEM from three, (C) from five, and (D) from four independent experiments. \*\*p≤0.01 and \*\*\*p≤0.001 compared to the incision activity of NEIL1 alone using one-way ANOVA and Tukey's post-hoc test. (C) NEIL1 was incubated with buffer or PARP-1 with or without NAD<sup>+</sup> and incubated with a non-radiolabeled oligonucleotide containing the 5OHU lesion as above. Samples were separated by SDS-PAGE and probed with anti-PAR, anti-PARP-1 and anti-NEIL1 antibodies.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Coordination of DNA repair by NEIL1 and PARP-1: a possible link to aging. *Aging (Albany NY)* (2012)



PARP-1 inhibits NEIL1 incision activity(A) NEIL1 (38 nM; 33 ng) was incubated with buffer, NAD<sup>+</sup> with or without PARP-1 (180 nM; 400 ng), or with PAR and reacted with a 5'- 32P-labeled oligonucleotide duplex containing a 5OHU lesion for 15 min at 37°C. The cleavage products were analyzed on a 20% denaturing gel containing 7 M urea. Percent incision was calculated by normalizing the amount of cleaved substrate (bottom band) to the amount of uncleaved product (top band). Data was then normalized to the amount of incision activity of NEIL1 alone (100%). (B) Incision assays were performed as in (A) in the presence of the indicated concentrations of PARP-1 with (activated PARP-1) or without NAD<sup>+</sup> (PARP-1) and quantified. (C) NEIL1 (33 ng) incision assays were performed in the presence of 400 ng of the indicated PARP-1 domains or full-length PARP-1 in the presence of NAD<sup>+</sup> (10 μM). (D) PARP-1 inhibits NEIL1 activity at both 5 nM and 20 nM concentrations of DNA substrate. The histograms in (A) and (B) represent the mean + SEM from three, (C) from five, and (D) from four independent experiments. \*\*p≤0.01 and \*\*\*p≤0.001 compared to the incision activity of NEIL1 alone using one-way ANOVA and Tukey's post-hoc test. (C) NEIL1 was incubated with buffer or PARP-1 with or without NAD<sup>+</sup> and incubated with a non-radiolabeled oligonucleotide containing the 5OHU lesion as above. Samples were separated by SDS-PAGE and probed with anti-PAR, anti-PARP-1 and anti-NEIL1 antibodies.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Coordination of DNA repair by NEIL1 and PARP-1: a possible link to aging. *Aging (Albany NY)* (2012)

## Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Poly(ADP-ribose) polymerase-1
Application	ELISA, ICC, WB
Application Notes	Detects bands ~116kDa (intact PARP) and ~85kDa (apoptosis-induced cleavage fragment) by Western blot.
Clone	C-2-10
Crossreactivity	Does not cross-react with chicken PARP.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Purified calf thymus poly(ADP-ribose) polymerase (PARP).
Isotype	IgG1
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1µg/ml, alkaline phosphatase)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from ascites.
Species Reactivity	Bovine, Hamster, Human, Monkey, Mouse, Rat



<b>Specificity</b>	Recognizes an epitope in the C-terminal part of the DNA binding domain of PARP.
<b>UniProt ID</b>	P18493
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)