

PARG (bovine thymus)

Citations: 11

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-202-0450	0.1U
--------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. For further dilution use the following enzyme dilution buffer: 50mM KPO ₄ , pH 7.5, 50mM KCl, 10 mM β-mercaptoethanol, 10% glycerol, 1mM DTT, 0.1% Triton.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Poly(ADP-ribose) glycohydrolase
Formulation	Liquid. In 50mM K-PO ₄ buffer, pH 7.5, containing 50mM potassium chloride, 0.1% Triton-x 100, 10% glycerol and 1mM DTT, 10mM β-Mercaptoethanol.
Purity Detail	Purified by substrate chromatography. The enzyme is composed of two major bands at 74kDa and 59kDa.
Source	Isolated from calf thymus.
Specific Activity	~20'000U/mg protein. 1U is defined as the amount of enzyme that hydrolyzes 1nmol of poly(ADP-ribose) per min. at 37°C.
UniProt ID	O02776

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)