

# 3-Aminobenzamide

## PARP inhibitor

Inhibits endogenous poly-ADP-ribosyltransferases often present in membrane preparations with minimal effect of bacterial toxin mediated ADP-ribosylation. Inhibits nitric oxide-induced apoptosis but not differentiation in HL-60 cells. Protects cells from oxygen radical and nitric oxide toxicity. Inhibits stress-induced apoptosis.

Citations: 12

[View Online »](#)

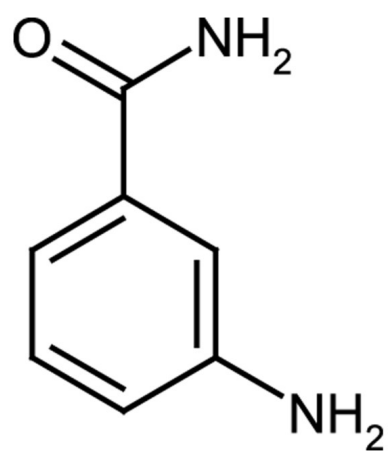
## Ordering Information

[Order Online »](#)

ALX-270-044-G001	1g
------------------	----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** Ambient

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 3-AB

**Appearance** White to tan powder.

**CAS** 3544-24-9

**Couple Target** PARP

**Couple Type** Inhibitor

**Formula**  $C_7H_8N_2O$

**MW** 136.2

**Purity**  $\geq 97\%$  (GC)

**RTECS** CU8992000

**Solubility** Soluble in DMSO or 100% ethanol.

**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)