

# Amifostine

HIF-1 $\alpha$  activator

Amifostine is a broad-spectrum radio- and chemoprotector. It restores transcriptional activity of specific p53 mutants, prevents chemotherapy-induced apoptosis and upregulates HIF-1 $\alpha$ .

Citations: 1

[View Online »](#)

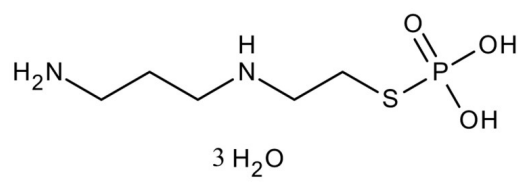
Ordering Information

[Order Online »](#)

|                |      |
|----------------|------|
| BML-GR341-0010 | 10mg |
|----------------|------|

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Long Term Storage** +4°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Appearance** White to off-white crystalline powder.

**CAS** 112901-68-5

**Couple Target** HIF

**Couple Type** Activator

**Formula**  $C_5H_{15}N_2O_3PS \cdot 3H_2O$

**Identity** Determined by IR.

**MW** 268.3

**Purity** ≥98% (UHPLC)

**Solubility** Soluble in water.

**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)