

# HBED

## Potent iron chelator

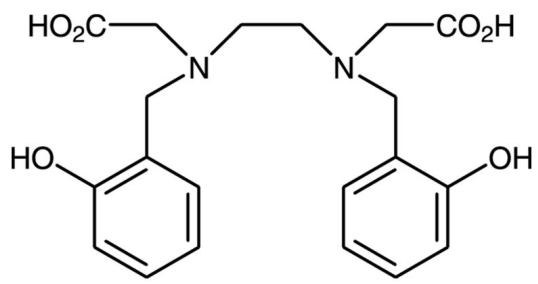
Iron chelator. Inhibits ferric ion catalyzed formation of hydroxyl radicals. A very potent ligand for the chelation of ferric iron *in vivo* (active in rat models but inactive in primate models). Inhibits neutrophil apoptosis. An important tool for probing the involvement of hydroxyl radicals in cellular processes.

Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



$\cdot\text{HCl}\cdot\text{H}_2\text{O}$

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store solutions at 20°C for up to 3 months.
<b>Long Term Storage</b>	Ambient
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Appearance</b>	White solid.
<b>CAS</b>	35369-53-0
<b>Formula</b>	$C_{20}H_{24}N_2O_6 \cdot HCl \cdot H_2O$
<b>MW</b>	442.9
<b>Purity</b>	≥95% (NMR)
<b>Solubility</b>	Soluble in DMSO (25mg/ml), in ethanol at 40°C (25mg/ml), in water (4mg/ml) at 50°C with vigorous agitation (approximately 15 minutes).

**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)