

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 »](#)

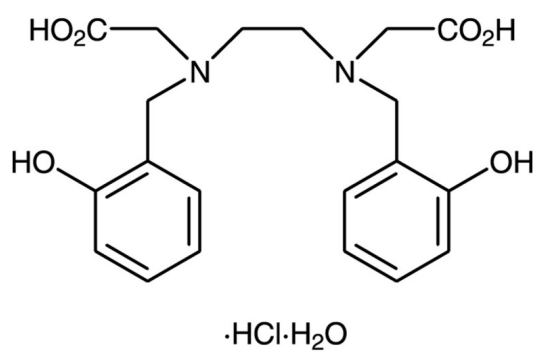
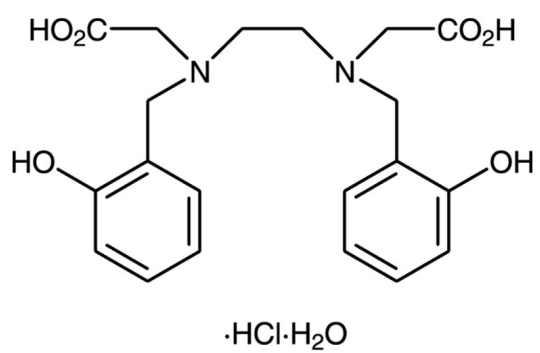
Ordering Information

[Order Online »](#)

BML-FR113-0250	250mg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	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.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	35369-53-0
Formula	$C_{20}H_{24}N_2O_6 \cdot HCl \cdot H_2O$
MW	442.9
Purity	≥95% (NMR)
Solubility	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)