

DTT

Protective agent for SH groups

Cell permeable protective agent for SH groups, maintaining monothiols completely in the reduced state and of reducing disulfides quantitatively. ER-stress inducer.

Citations: 14

[View Online »](#)

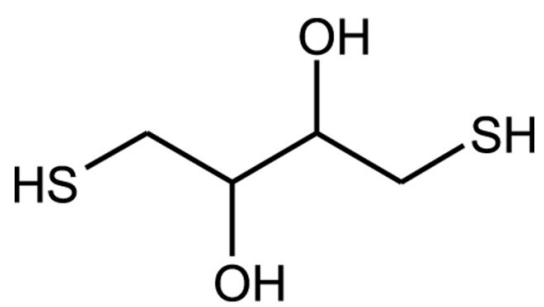
Ordering Information

[Order Online »](#)

ALX-280-001-G005	5g
ALX-280-001-G010	10g
ALX-280-001-G025	25g
ALX-280-001-G050	50g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Very hygroscopic. Keep under inert gas.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1,4-Dithio-DL-threitol, DL-Dithiothreitol, Cleland's Reagent
Appearance	White to off-white solid.
CAS	3483-12-3
Couple Type	Ligand
Formula	$C_4H_{10}O_2S_2$
Identity	Determined by IR.
MI	14: 3376
MW	154.3
Purity	≥98% (Assay by Iodometric Titration)
RTECS	EK1610000
Solubility	Soluble in water, 100% ethanol, acetone or ethyl acetate.

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-
fr@enzolifesciences.com](mailto:info-fr@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)