

2-

Mercaptoethanesulfonate Sodium

Organosulfur, antioxidant.

A uroprotective agent used in combination with cancer chemotherapeutic drugs to decrease urotoxicity. It was found to have antitumor activity against several cell lines. In addition, it is used as an antimucolytic agent.

Ordering Information

[Order Online »](#)

LKT-M1774-G025	25g
----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2-Mercaptoethanesulfonic acid sodium salt, MESNA

CAS 19767-45-4

Formula $C_2H_5O_3S_2 \cdot Na$

MI 13: 5936

MW 164.18

Purity $\geq 98\%$

Solubility Soluble in water (50 mg/mL), DMSO (33 mg/mL). Slightly soluble in alcohol. Practically insoluble in cyclohexane.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com