

1,3,4,5-Tetramethyl-2- [(2-oxo- propyl)thio]imidazolium . chloride

1,3,4,5-Tetramethyl-2-[(2-oxopropyl)thio]imidazolium . chloride is a potent, active site directed inhibitor of transglutaminases.

Ordering Information

[Order Online »](#)

NZY-T101-010	10mg
--------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name L-682,777, 1,3,4,5-Tetramethyl-2-(2-oxo-propylsulfanyl)-3H-imidazol-1-ium . Cl

Appearance Off-white solid.

Formula $C_{10}H_{17}N_2OS \cdot Cl$

MW 213.3 . 35.5

Purity ≥95% (HPLC)

Reconstitution Predissolve in DMSO, then dilute with aqueous buffers.

Solubility Soluble in DMSO, dimethyl formamide; poorly soluble in methanol.

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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