

THPTA (25 mg)

Ordering Information

[Order Online »](#)

JBS-CLK-1010-25	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tris((1-hydroxy-propyl-1H-1,2,3-triazol-4-yl)methyl)amine

Appearance off-white to grey solid

CAS 760952-88-3

Formula $C_{18}H_{30}N_{10}O_3$

MW 434.5

Purity > 95 % (HPLC)

Solubility water, DMSO, DMF, MeOH

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