

Guanosine 5'-O-(3-thiotriphosphate) . tetralithium salt

Non-hydrolyzable GTP analog

Guanosine 5'-O-(3-thiotriphosphate) (GTP-g-S) is a non-hydrolyzable analog of guanosine 5'-O-triphosphate (GTP) (Prod. No. ALX-480-057). GTP-g-S stimulates inositol trisphosphate formation in membranes.

Citations: 3

[View Online »](#)

Ordering Information

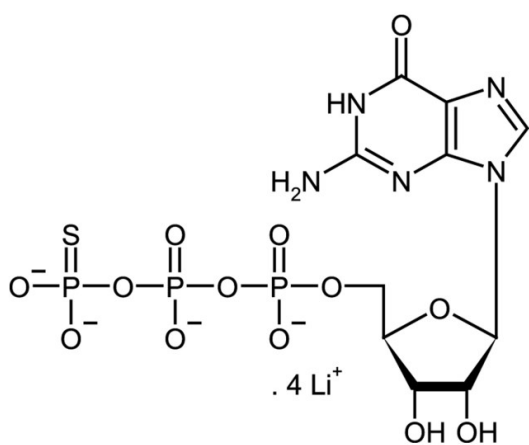
[Order Online »](#)

ALX-480-056-M001	1mg
ALX-480-056-M005	5mg

Manuals, SDS & CofA

[View Online »](#)

- Used to obtain crystal structures of GTPases and measure their stability in presence of ligands



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name GTP-γ-S . 4Li

Appearance White to off-white solid.

CAS 94825-44-2

Formula $C_{10}H_{12}N_5O_{13}P_3S \cdot 4Li$

MW 535.2 . 27.8

Purity ≥85% (HPLC)

Solubility Soluble in water (25mg/ml).

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)