

Biotin-isoxazole

Precipitates RNA Binding Proteins

Exposure of cell or tissue lysates to biotin-isoxazole precipitates hundreds of RNA-binding proteins with significant overlap to the constituents of RNA granules. In one study roughly 580 nuclear proteins were identified as being precipitated by B-Isox including TAF15, RNA polymerase II (largest subunit), numerous subunits of the mediator complex as well as a variety of enzymes involved in epigenetic modification of histones. SR domains were found to be the determinant that facilitates B-Isox precipitation. Also allows for the fractionation / isolation of SG-associated proteins and RNA precipitates.

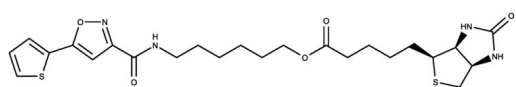
Ordering Information

[Order Online »](#)

ENZ-CHM500-0005	5mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



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 6-(5-(Thiophen-2-yl)isoxazole-3-carboxamido)hexyl 5-((3aS,4S,6aR)-2-oxohexahydro-1H-thieno[3,4-d]imidazol-4-yl)pentanoate

Appearance Beige solid.

CAS 1377605-22-5

Formula $C_{24}H_{32}N_4O_5S_2$

Identity Determined by NMR.

MW 520.66

Purity ≥98% (TLC)

Solubility Soluble in DMSO (30 mg/ml).



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)