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

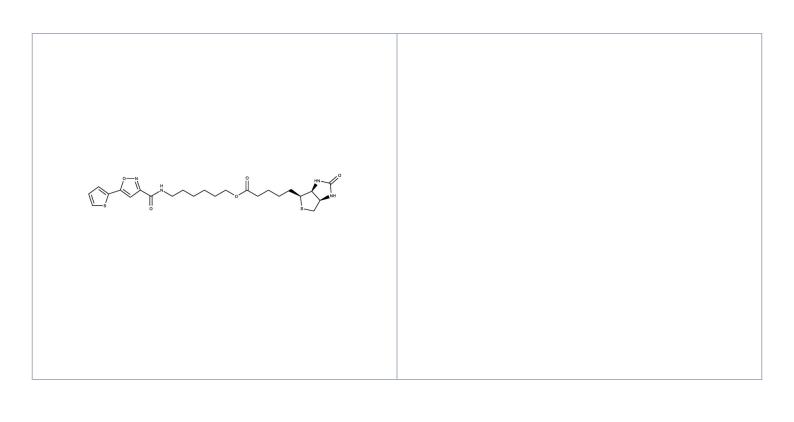
Order Online »

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).

Last modified: February 27, 2025