

dBET6

Optimized and highly cell-permeable
chemical degrader of BET bromodomain
proteins, based on PROTAC

dBET6 is chemical degrader of BET bromodomain proteins based on PROTAC and led a collapse of global elongation that phenocopies CDK9 inhibition.

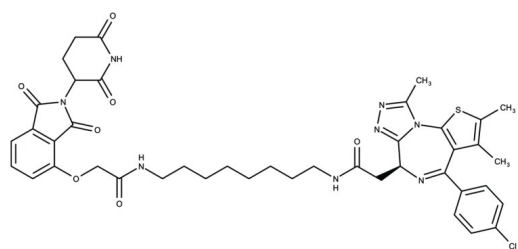
Ordering Information

[Order Online »](#)

ENZ-CHM257-0001	1mg
ENZ-CHM257-0005	5mg
ENZ-CHM257-0010	10mg
ENZ-CHM257-0050	50mg

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to light yellow solid.

CAS 1950634-92-0

Couple Target BET bromodomain protein, Ligase [E3]

Couple Type Inhibitor, Ligand

Formula $C_{42}H_{45}ClN_8O_7S$

MW 841.4

Purity ≥98% (LCMS)

Solubility Soluble in DMSO (240 mg/mL). Insoluble in water.

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-
be@enzolifesciences.com](mailto:info-be@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)