

A1874

Nutlin-based PROTAC targeting on BRD4

A1874 is a nutlin-based PROTAC targeting on BRD4, consist of a BRD4-ligand JQ1 linked to a MDM2 E3 ligase ligand RG7388. A1874 possesses dual function as it dose-dependently degraded its target BRD4, accompanied with the downregulation of c-MYC, as well as stabilization of P53 and upregulated P21 due to the efficient engagement of MDM2 by RG7388.

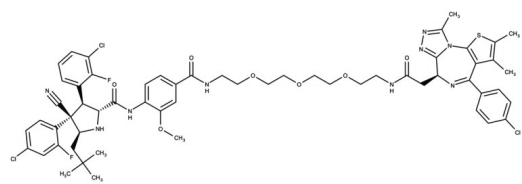
Ordering Information

[Order Online »](#)

ENZ-CHM262-0005	5mg
ENZ-CHM262-0010	10mg
ENZ-CHM262-0025	25mg

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 solid.
CAS	2064292-12-0
Couple Target	BET bromodomain protein, Ligase [E3]
Couple Type	Inhibitor, Ligand
Formula	$C_{58}H_{62}Cl_3F_2N_9O_7S$
MW	1173.6
Purity	≥98%

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)