

Arv-825

PROTAC targeting on BRD4

ARV-825 is a PROTAC targeting on BRD4, consisting of a BRD4-ligand JQ-1 linked to a cereblon E3 ligase ligand Pomalidomide, with K_d s values of 90nM and 28nM for affinity to BD1 and BD2.

Ordering Information

[Order Online »](#)

ENZ-CHM258-0001	1mg
-----------------	-----

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	Yellow solid.
CAS	1818885-28-7
Couple Target	BET bromodomain protein, Ligase [E3]
Couple Type	Inhibitor, Ligand
Formula	$C_{46}H_{47}ClN_8O_9S$
MW	923.4
Purity	≥98%

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com