

Compound 115-7c

HSP70 modulator

Acts as an artificial co-chaperone for HSP70. The activity of compound 115-7c mirrors those of HSP40 stimulating the ATPase and protein-folding activities of prokaryotic HSP70. Increases tau levels in cell lines overexpressing tau. Induces solubilization of polyglutamine and exacerbates polyQ-mediated toxicity. An important probe for understanding HSP70 functions.

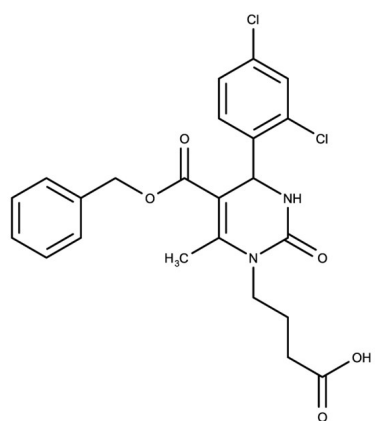
Ordering Information

[Order Online »](#)

ENZ-CHM471-0025	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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 4-(5-((Benzyloxycarbonyl)-4-(2,4-dichlorophenyl)-6-methyl-2-oxo-3,4-dihydropyrimidin-1(2H)-yl)butanoic acid

Appearance Off-white solid.

CAS 908074-72-6

Couple Target HSP70

Couple Type Activator

Formula $C_{23}H_{22}Cl_2N_2O_5$

Identity Determined by NMR.

MW 477.34

Purity ≥98% (TLC)

Solubility Soluble in DMSO (20 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)