

# 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.

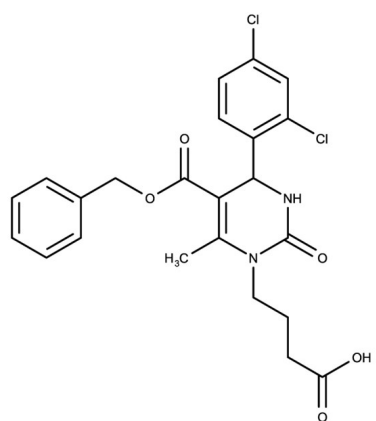
### Ordering Information

[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).

**Last modified: February 27, 2025**



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-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@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)