

SMER28

**Small molecule inducer of autophagy
important for neurodegenerative disease
research.**

Chemical inducer of autophagy which acts via an mTOR-independent mechanism. It increased autophagosome synthesis and enhanced the clearance of model autophagy substrates such as A53T α -synuclein and mutant huntingtin fragments. It reduced A β peptide and AAP-CTF levels in an Alzheimers model and attenuated mutant huntingtin-fragment toxicity in Huntington's disease models, suggesting therapeutic potential.

Citations: 4

[View Online »](#)

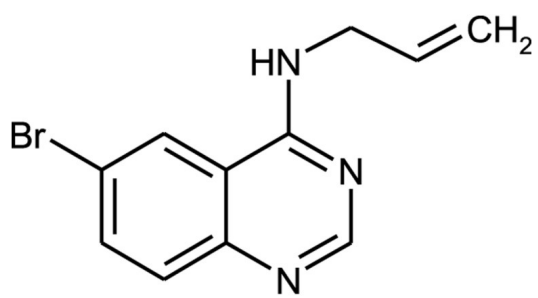
Ordering Information

[Order Online »](#)

BML-EI397-0005	5mg
BML-EI397-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	6-Bromo-4-allylaminoquinazoline
Appearance	Off-white solid.
CAS	307538-42-7
Formula	$C_{11}H_{10}BrN_3$
MW	264.1
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (8mg/ml) or 100% ethanol (2mg/ml).
Source	Synthetic.



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