

Thiazovivin

ROCK inhibitor

Thiazovivin is a potent and selective inhibitor of Rho-associated kinase (ROCK), an enzyme involved in regulating cell shape, motility, and differentiation. It is widely used in cell culture techniques, particularly for the generation and maintenance of induced pluripotent stem cells (iPSCs).

Ordering Information

[Order Online »](#)

ENZ-CHM322-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

- GMP format available

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Powder: Stable at -20°C for 3 years or +4°C for 2 years; In solvent: Stable at -80°C for 2 years or -20°C for 1 year.
Handling	Before opening the vial, please centrifuge to gather the compound at the bottom of the vial.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	N-(phenylmethyl)-2-(4-pyrimidinylamino)-4-thiazolecarboxamide, N-Benzyl-2-(pyrimidin-4-ylamino)thiazole-4-carboxamide
Appearance	Light yellow to brown solid.
CAS	1226056-71-8
Couple Target	ROCK
Couple Type	Inhibitor
Formula	$C_{15}H_{13}N_5OS$
Identity	Identity determined by LCMS.
MW	311.36
Purity	≥99% (LCMS)
Solubility	Soluble in DMSO (30 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-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)