

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

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

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

Last modified: November 7, 2024