

Product Information

Description

ROCK inhibitor. Increases the survival of cells differentiated from induced pluripotent stem cells upon thawing. Greatly enhances the efficiency of fibroblast reprogramming to generate induced pluripotent stem cells (iPSCs) when used in combination with SB 431542 and PD 0325901. Improves the survival of human embryonic stem cells (hESCs) and human induced pluripotent stem cells (hiPSCs) upon dissociation.

Target:

ROCK

Chemical Name:

N-benzyl-2-(pyrimidin-4-ylamino)-1,3-thiazole-4-carboxamide

Molecular Formula:C₁₅H₁₃N₅OS**Supplied As:**

Solid

Solubility:

63 mg/mL in DMSO, insoluble in water, insoluble in ethanol.

Stability:

At least 6 months at -20°C. Anhydrous DMSO stock solutions are stable for at least 1 year at -80°C and 1 month at -20°C. Avoid freeze/thaw cycles.

Storage:

-20°C

MW:

311.36 Da

Cas No.

1226056-71-8

Purity:

≥99%

Chemical Structure

