

Aurora A Inhibitor I

(Catalog No. M75000)

Alternative Names

4-(5-Amino-1-(2,6-difluorobenzoyl)-1H-[1,2,4]triazol-3-ylamino)-benzenesulfonamide

Description

4-(5-Amino-1-(2,6-difluorobenzoyl)-1H-[1,2,4]triazol-3-ylamino)-benzenesulfonamide is a cell-permeable triazolylsulfonamido compound that acts a reversible, ATP-competitive kinase inhibitor with selectivities towards Aurora kinases (IC₅₀ = 11 and 15 nM for Aurora-A, Aurora-B, respectively) and Cdk's (IC₅₀ = 9, 4, and 3 nM for Cdk1/B, Cdk2/A, and Cdk2/E, respectively). Displays antitumor properties both in vitro (IC₅₀ in the range of 112 - 514 nM against various human cancer cell lines) and in a xenograft murine model in vivo (75-100 mg/kg, i.p.).

CAS Number

443797-96-4

Chemical Formula

C₁₅H₁₂F₂N₆O₃S • CH₃CN

Molecular Weight

435.4

Size

5 mg

Purity

≥97%

Solubility

Soluble in DMSO

Storage Temp

-20°C

Ordering Information:

Products

Aurora A Inhibitor I

Size

5 mg

Cat. No.

M75000 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.