

HA-1077

Catalog #: 27030

Lot #: 140418

Size: 250 mg

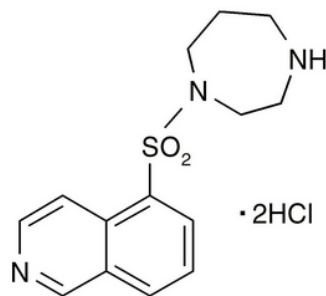
Structure:

CAS Registry #: 203911-27-7

Purity: ≥99%

Chemical Formula: C₁₄H₁₇N₃O₂S·2HCl

Molecular Weight: 364.3



Description: Dihydrochloride form of HA-1077. Inhibits a number of kinases, including ROCK I (IC₅₀ = 10.7 μM), ROCK II (IC₅₀ = 1.9 μM), MAPK1 (IC₅₀ = 5 μM), and PKCRK2 (IC₅₀ = 4 μM). Weaker inhibitor of PKC (IC₅₀ = 425 μM) and MLCK (IC₅₀ = 95 μM).

Synonym(s): Fasudil, Dihydrochloride Salt, 1-(5-Isoquinolinesulfonyl)homopiperazine, 2HCl, Eril, AT-877

Appearance: White crystalline powder.

Solubility: Very soluble in water, up to 200 mg/ml or more; soluble in DMSO up to about 10 mg/ml; not soluble in ethanol; addition of an aqueous solution of this dihydrochloride salt into a buffer solution may significantly alter its solubility

Storage/Stability: Store at or below 20°C. Solid form is stable at least 12 months from date of receipt, when stored as directed. Do not store aqueous solutions for more than one day.

References:

1. Davies, S.P., *et al.*, *Biochem. J.* 2000;**351**:95-105.