HA-1077

Rho kinase (ROCK) inhibitor

Inhibits Rho kinase (ROCK; K_i =0.33 μ M). Selective protein kinase A and G inhibitor (both K_i 's = 1.6 μ M). Inhibits agonist-induced activation of NADPH oxidase in human neutrophils (IC $_{50}$ =15 μ M). Inhibitor of myosin light chain kinase and Ca $^{2+}$ /calmodulin-dependent protein kinase II. Inhibits translocation of PKC β I, PKC β II and PKC ζ . Cell permeable Ca $^{2+}$ antagonist with antivasospastic properties.

Citations: 14

View Online »

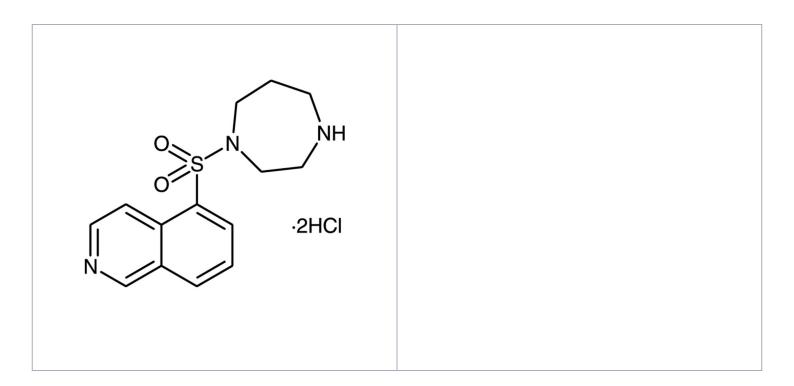
Ordering Information

Order Online »

BML-EI233-0010	10mg
BML-EI233-0050	50mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Stable for at least

1 year after receipt when stored, as supplied, at 0-4°C. Stock solutions are stable for up

to 3 months at -20°C.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1-(5-Isoquinolinesulfonyl)homopiperazine . 2HCl, Fasudil

Appearance White to off-white solid.

CAS 103745-39-7

Couple Target CaM kinase, MLCK, PKA, PKG, ROCK

Couple Type Inhibitor

Formula $C_{14}H_{17}N_3O_2S$. 2HCl

Identity Determined by 1H-NMR and MS.

MI 14: 3942

MW 364.3

MeltingPoint 238-242°C

Purity ≥98% (TLC)

Solubility Soluble in water (>20mg/ml), methanol or DMSO.