

Product Information

Description

Allosteric and reversible inhibitor of protease-activated receptor 1 (PAR1)-mediated platelet activation ($IC_{50} = 0.26 \mu M$ for the inhibition of platelet P-selectin expression on human platelets).

Target:

PAR1

Pathway

Angiogenesis

Chemical Name:

2-bromo-N-[3-(butanoylamino) phenyl]benzamide

Molecular Formula:

$C_{17}H_{17}BrN_2O_2$

Supplied As:

Solid

Solubility:

insoluble in H_2O ; ≥ 16.4 mg/ml in DMSO; ≥ 46.9 mg/ml in EtOH with ultrasonic. For obtaining a higher solubility, please warm the tube at $37^\circ C$ and shake it in the ultrasonic bath for a while. Long-term storage in solution is not recommended. Not stable in diluted solutions.

Stability:

At least 6 months at $-20^\circ C$. Avoid freeze/thaw cycles.

Storage:

$-20^\circ C$

MW:

361.23 Da

Cas No.

423735-93-7

Purity:

$\geq 97\%$

Chemical Structure

