PF-622

FAAH inhibitor

Novel, potent and selective inhibitor of fatty acid amide hydrolase (FAAH). Inhibits FAAH in a time-dependent manner by covalently modifying the enzyme's active site serine (IC_{50} =33nM). It is completely selective for FAAH relative to other mammalian serine hydrolases.

Citations: 1

View Online »

Ordering Information

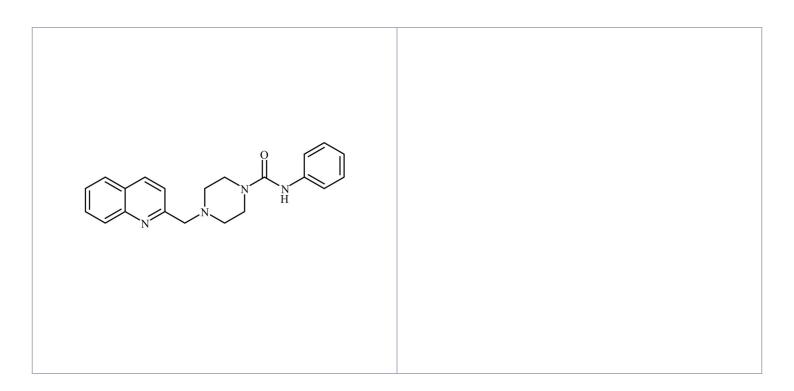
Order Online »

BML-EI402-0005

5mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Store, as supplied,

at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-phenyl-4-(quinolin-2-ylmethyl)piperazine-1-carboxamide

Appearance Off-white solid.

CAS 898235-65-9

Couple Target FAAH

Couple Type Inhibitor

Formula $C_{21}H_{22}N_4O$

MW 346.4

Purity ≥98% (TLC)

Solubility Soluble in DMSO (7mg/ml) or 100% ethanol (3mg/ml).

Last modified: May 29, 2024