URB-597

FAAH inhibitor

URB-597 is a potent and selective inhibitor of the enzyme fatty acid amide hydrolase (FAAH), which is involved in the hydrolysis of endogenous cannabinoids such as an anadamide. Inhibition of FAAH may serve to enhance the analgesic effects of endogenous cannabinoids. URB-597 has been shown to inhibit FAAH in rat cortical neuron membranes and in intact neurons with IC $_{50}$ values of 4.6nM and 0.5nM, respectively. The compound has also been shown to have anxiolytic effects in rats (ID $_{50}$ =0.15mg/kg). Because URB-597 acts on FAAH and does not interfere with the binding of an anadamide to the CB1 and CB2 receptors, it is a novel tool for drug development.

Citations: 6

View Online »

Ordering Information

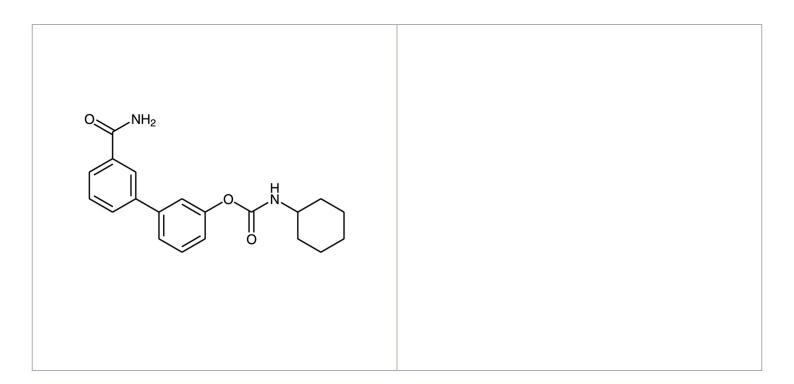
Order Online »

BML-EI327-0025

25mg

Manuals, SDS & CofA

View Online »



Handling & Storage

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

at -20°C for up to 1 year. Store solutons at -20°C for up to 3 months.

Handling Protect from light.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Cyclohexylcarbamic acid 3'-carbamoyl-biphenyl-3-yl ester

Appearance White solid.

CAS 546141-08-6

Couple Target FAAH

Couple Type Inhibitor

Formula $C_{20}H_{22}N_2O_3$

MW 338.4

MeltingPoint 155-157°C

Purity ≥98% (TLC)

Solubility Soluble in 100% ethanol (1mg/ml), dimethyl formamide

(10mg/ml) or DMSO (10mg/ml).

Last modified: May 29, 2024



uk@enzolifesciences.com