# **URB-602**

## Monoacylglycerol lipase inhibitor

Selectively inhibits monoacylglycerol (MAG) lipase (IC $_{50}$ =28 $\mu$ M) via a noncompetitive mechanism. Does not inhibit FAAH, DAG lipase or COX-2. URB-602 selectively raises levels of 2-arachidonoylglycerol with no change in anandamide concentrationsin cells and tissues. In the same way anandamide levels were markely elevated by the FAAH inhibitor URB597 at 1 $\mu$ M.

Citations: 1

View Online »

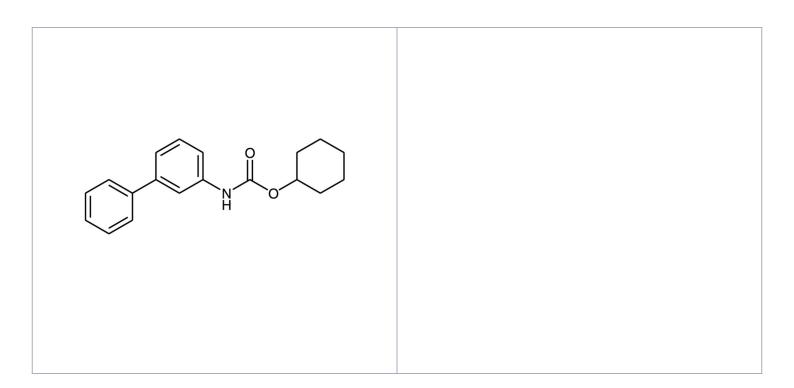
**Ordering Information** 

Order Online »

**BML-EI348-0050** 50mg

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 Biphenyl-3-yl carbamic acid cyclohexyl ester

**Appearance** Off-white solid.

**CAS** 565460-15-3

Couple Target Monoacylglycerol lipase

Couple Type Inhibitor

Formula  $C_{19}H_{21}NO_2$ 

**MW** 295.4

MeltingPoint 122-123°C

Purity ≥98% (TLC)

**Solubility** Soluble in DMSO (30 mg/ml) or ethanol (5 mg/ml warm).

Last modified: May 29, 2024

