

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 anandamide. 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 anandamide to the CB1 and CB2 receptors, it is a novel tool for drug development.

Citations: 6

[View Online »](#)

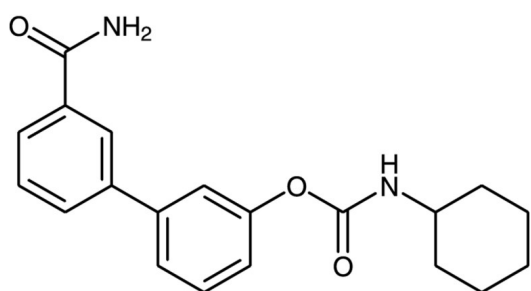
Ordering Information

[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 solutions 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).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com