

U-37883A

K⁺ channel blocker

A non-sulfonylurea inhibitor of K-ATP channels. Inhibits vascular K-ATP channels. Inhibits dopamine-modulated potassium channels.

Citations: 2

[View Online »](#)

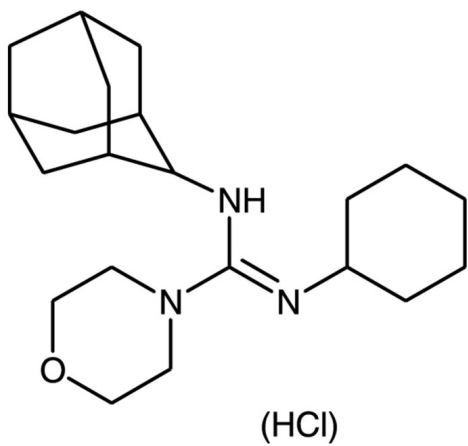
Ordering Information

[Order Online »](#)

BML-KC159-0005	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 1 month.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	N-(1-Adamantyl)-N'-cyclohexyl-4-morpholinecarboxamide . HCl
Appearance	White to off-white solid
CAS	57568-80-6
Couple Target	Potassium channel
Couple Type	Blocker
Formula	$C_{21}H_{35}N_3O \cdot HCl$
MW	382
Purity	≥98% (TLC)
Solubility	Soluble in water (5mg/ml).

Last modified: May 29, 2024



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-be@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