

Acetylcholine receptor activator

Highly selective allosteric potentiator of the M₁ muscarinic acetylcholine receptor (mAChR). Reduces the concentration of ACh required to activate M₁ up to 129-fold and displays no potentiation, agonism or antagonism activity at other mAChRs at concentrations up to 100μM. Reverses scopolamine-induced memory deficits and increases blood flow to the cerebral cortex in mouse models. M₁ activators slow the progression of Alzheimers disease. Induces β-arrestin recruitment to M₁. At 100μM, it activated M₁ in the absence of ACh to ca. 50% maximal response. At 10μM, it reduced the concentration of ACh required to displace radioligand by 45-fold. *In vivo* experiments were carried out in mice at 15 or 20mg/kg. The compound is cell permeable, brain penetrant and orally active.

Citations: 3

[View Online »](#)

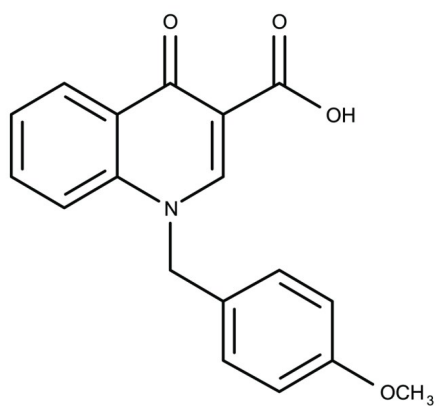
Ordering Information

[Order Online »](#)

BML-C121-0010	10mg
BML-C121-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at room temperature. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-(4-Methoxybenzyl)-4-oxo-1,4-dihydro-3-quinoline carboxylic acid, Benzyl quinolone carboxylic acid
Appearance	White solid.
CAS	338747-41-4
Couple Target	Acetylcholine receptor
Couple Type	Activator
Formula	$C_{18}H_{15}NO_4$
MW	309.3
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO (10mg/ml).
Source	Synthetic.



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