

Icilin

TRPM8 activator

Cooling agent. Strongly activates TRPM8 ($IC_{50}=0.2\mu M$), cold menthol receptor 1 (CMR1, $IC_{50}=0.36\mu M$), a nonselective, cold-activated cation channel and TRPA1 (at 10- to 100-fold higher concentration). Induces currents in TRPM8 expressing HEK 293 cells ($EC_{50}=0.36mM$) more potently than menthol (Prod. No. ALX-420-042) or low temperatures. Activates TRPM8 with variable latency, followed by extensive desensitization provided calcium is present. The effects of icilin can be impacted by both pH and cytosolic calcium levels.

Citations: 9

[View Online »](#)

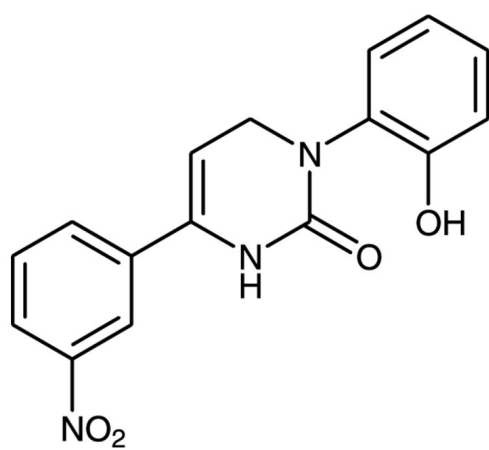
Ordering Information

[Order Online »](#)

BML-CM126-0050	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	AG 3-5, 3,6-Dihydro-1-(2-hydroxyphenyl)-4-(3-nitrophenyl)-2(1H)-pyrimidinone
Appearance	Light yellow to orange solid.
CAS	36945-98-9
Couple Target	TRP channel
Couple Type	Activator
Formula	$C_{16}H_{13}N_3O_4$
MW	311.3
Purity	≥97% (TLC)
Solubility	Soluble in DMSO (50mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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