

Adapalene

Retinoic acid receptor ligand

Adapalene is a stable analog of retinoic acid and acts selectively as an agonist at RAR β and RAR γ . It inhibits proliferation and induces apoptosis in colorectal cancer cells. Its unique pharmacological properties make it superior to other retinoids for the treatment of acne.

Citations: 3

[View Online »](#)

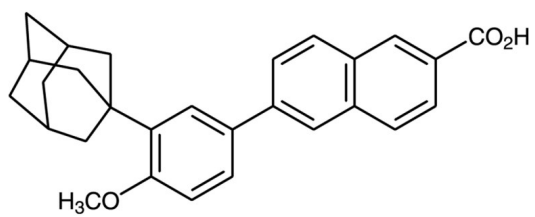
Ordering Information

[Order Online »](#)

BML-GR111-0020	20mg
----------------	------

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	6-[3-(1-Adamantyl)-4-methoxyphenyl]-2-naphthoic acid
Appearance	White solid.
CAS	106685-40-9
Couple Target	Retinoic acid receptor
Couple Type	Ligand
Formula	$C_{19}H_{34}O_3$
MW	412.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (3 mg/mL) or DMF (6 mg/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-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