

5-Ketoeicosa- 6E,8Z,11Z,14E- tetraenoic acid

R527 ligand

Stimulates neutrophil degranulation and Ca^{2+} mobilization by a mechanism independent of the LTB₄ receptor. Has been shown to act as an agonist at a G protein-coupled receptor (R527). Induces the infiltration of granulocytes into human skin.

Citations: 6

[View Online »](#)

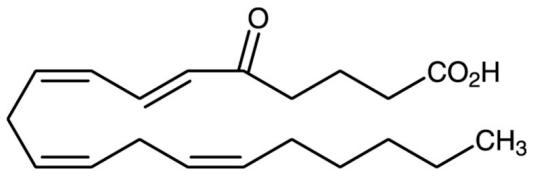
Ordering Information

[Order Online »](#)

BML-K005-0050	50µg
---------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -80°C for up to 1 year.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Liquid.
CAS	106154-18-1
Couple Target	R527
Couple Type	Ligand
Formula	$C_{20}H_{30}O_3$
Formulation	Oil dissolved in ethanol
MW	318.5
Purity	≥95% (HPLC)

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