

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

## R527 ligand

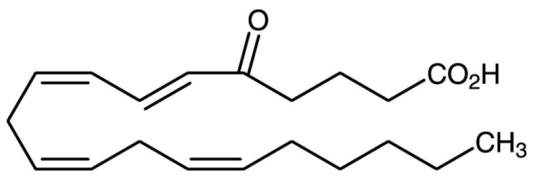
Stimulates neutrophil degranulation and  $\text{Ca}^{2+}$  mobilization by a mechanism independent of the LTB<sub>4</sub> receptor. Has been shown to act as an agonist at a G protein-coupled receptor (R527). Induces the infiltration of granulocytes into human skin.

Citations: 6

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

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

**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-  
be@enzolifesciences.com](mailto:info-be@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)