

# MK-886

## FLAP inhibitor

Inhibitor of PPAR $\alpha$  activation. Inhibits 5-lipoxygenase-activating protein (FLAP) and leukotriene biosynthesis. Binds to FLAP with high affinity and prevents the activation of 5-lipoxygenase. IC<sub>50</sub>'s for leukotriene biosynthesis: 2.5 nM in intact leukocytes; 1.1  $\mu$ M in whole blood. Induces apoptosis through the mitochondrial pathway.

Citations: 9

[View Online »](#)

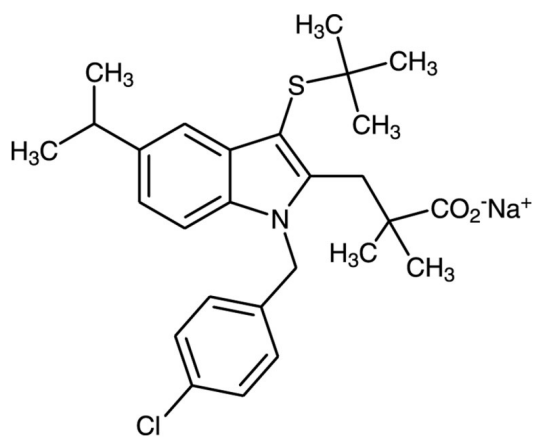
## Ordering Information

[Order Online »](#)

BML-EI266-0005	5mg
----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at room temperature. Prepare fresh solutions daily.
<b>Long Term Storage</b>	Ambient
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	L-663,536, 3-[1-(p-Chlorobenzyl)-5-(isopropyl)-3-t-butylthioindol-2-yl]-2,2-dimethylpropanoic acid . Na
<b>Appearance</b>	White solid.
<b>CAS</b>	118414-82-7
<b>Couple Target</b>	FLAP, PPAR
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{27}H_{33}ClNO_2S \cdot Na$
<b>MW</b>	494.0
<b>Purity</b>	≥98% (HPLC)
<b>Solubility</b>	Soluble in 100% ethanol (25mg/ml) or DMSO (25mg/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-  
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)