

LY 171883

Leukotriene D₄ antagonist

Selective, orally active leukotriene D₄ antagonist. Has been shown to bind to the γ -isoform of the peroxisome proliferator-activated receptor (PPAR γ).

Citations: 6

[View Online »](#)

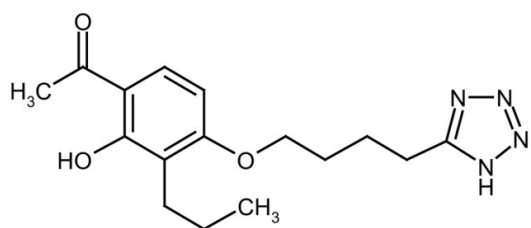
Ordering Information

[Order Online »](#)

BML-RA101-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Organic stock solutions are stable for at least six months when stored at -20°C.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tomelukast

Appearance White solid.

CAS 88107-10-2

Couple Target Leukotriene, PPAR, Prostanoid receptor

Couple Type Inhibitor, Ligand

Formula $C_{16}H_{22}N_4O_3$

MW 318.4

MeltingPoint 117-119°C

Purity ≥90% (HPLC)

Solubility Soluble in DMSO (25mg/ml), 100% ethanol (25mg/ml) or dimethyl formamide. Also soluble in 0.5M aqueous sodium bicarbonate.

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: +33 472 440 656
[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)