

MK-571 . sodium salt

Leukotriene receptor ligand

A potent, selective, orally active leukotriene D₄ antagonist. K_i=2.1 nM for inhibition of [³H]-LTD₄ binding to human lung membranes. Competitively antagonizes contractions of guinea pig trachea and ileum and human trachea induced by LTD₄. Blocks LTD₄ and ascaris-induced bronchoconstriction in conscious squirrel monkeys.

Citations: 31

[View Online »](#)

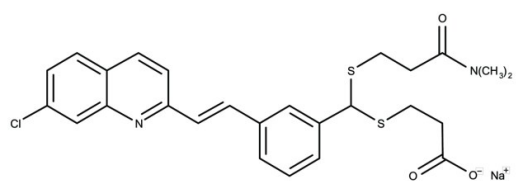
Ordering Information

[Order Online »](#)

BML-RA109-0005	5mg
BML-RA109-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 2 months at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	L660,711
Appearance	Off-white solid.
CAS	115104-28-4
Couple Target	Leukotriene receptor, Prostanoid receptor
Couple Type	Ligand
Formula	$C_{26}H_{26}ClN_2O_3S_2 \cdot Na$
MW	537
Purity	≥90% (HPLC)
Solubility	Soluble in water (20mg/ml), aqueous buffers, 100% ethanol (10mg/ml) or DMSO.

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: +33 472 440 655
[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)