

Tranilast

Anti-inflammatory and anti-angiogenic

Anti-inflammatory and anti-angiogenic agent. Inhibits VEGF-induced angiogenesis *in vivo*. Inhibits proliferation and tube formation of human endothelial cells *in vitro*. Also inhibits smooth muscle cell proliferation and prevents post-angioplasty restenosis. Synthetic compound with structural similarity to the tryptophan (Trp) catabolites kynurenines. Tranilast ameliorates the disease severity in animal models of multiple sclerosis, in part, by inducing T cell tolerance. Trp degradation is an important mechanism in the maintenance of T cell tolerance in a variety of disease models and is mediated by the rate limiting enzyme indoleamine 2,3-dioxygenase (IDO).

Citations: 11

[View Online »](#)

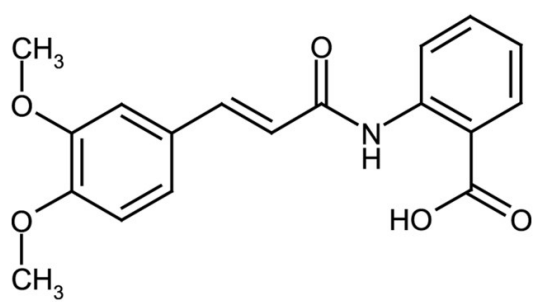
Ordering Information

[Order Online »](#)

ALX-550-409-M010	10mg
------------------	------

Manuals, SDS & CoFA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(3',4'-Dimethoxycinnamoyl)anthranilic acid, 3,4-DAA
Appearance	Off-white to yellow solid.
CAS	53902-12-8
Formula	$C_{18}H_{17}NO_5$
MI	14: 9570
MW	327.3
Purity	≥95% (HPLC)
RTECS	DG8731000
Solubility	Soluble in 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)