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-angiplasty retsenosis. 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

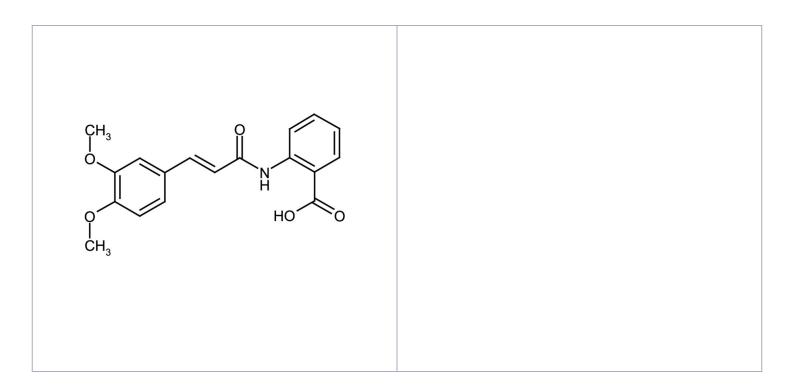
Order Online »

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