

Leflunomide

Immunosuppressant

Immunosuppressive drug that inhibits T and B cell responses *in vivo*. Inhibitor of tyrosine phosphorylation (tyrosine kinase inhibitor), of dihydroorotate dehydrogenase, an enzyme for *de novo* pyrimidine synthesis, of TNF-dependent NF- κ B activation, and of COX-2. The physiological activity of leflunomide is attributed mainly to its metabolite A77 1726 (Prod. No. ALX-430-096).

Citations: 22

[View Online »](#)

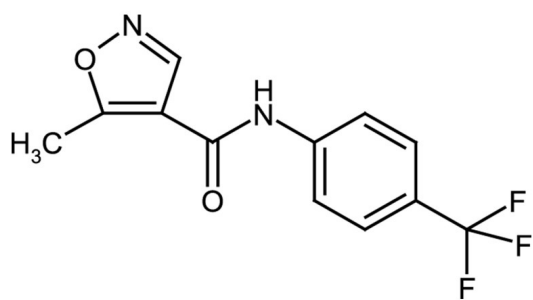
Ordering Information

[Order Online »](#)

ALX-430-095-M050	50mg
ALX-430-095-G001	1g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Keep under inert gas.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	HWA486, 5-Methylisoxazole-4-(4-trifluoromethylcarboxanilide), N-(4-Trifluoromethylphenyl)-5-methylisoxazol-4-carboxamide
Appearance	White to off-white crystalline powder.
CAS	75706-12-6
Couple Target	COX, NF-kappaB
Couple Type	Inhibitor
Formula	$C_{12}H_9F_3N_2O_2$
Identity	Determined by IR.
MI	14: 5432
MW	270.2
Purity	≥98.5% (HPLC)
RTECS	NY2354200
Solubility	Soluble in DMSO or methanol.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com