

# A77 1726

## Dihydrorotate dehydrogenase inhibitor

Physiologically active metabolite of the immunosuppressive drug leflunomide (Prod. No. ALX-430-095). Inhibits the activity of dihydrorotate dehydrogenase and of protein tyrosine kinases. Blocks TNF-mediated NF- $\kappa$ B activation in a dose- and time-dependent manner. Also inhibits the activity of cyclooxygenase-2 (COX-2) *in vitro* and *in vivo*.

Citations: 10

[View Online »](#)

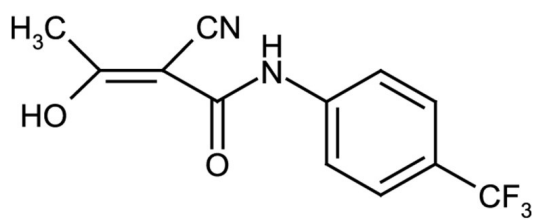
## Ordering Information

[Order Online »](#)

ALX-430-096-M005	5mg
ALX-430-096-M025	25mg

## 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	N-(4-Trifluoromethylphenyl)-2-cyano-3-hydroxycrotonamide, 2-Cyano-3-hydroxy-N-[4-(trifluoromethyl)phenyl]-2-butenamide
Appearance	White solid.
CAS	108605-62-5
Couple Target	Dihydrorotate dehydrogenase
Couple Type	Inhibitor
Formula	$C_{12}H_9F_3N_2O_2$
Identity	Determined by NMR.
MW	270.2
Purity	≥98% (HPLC)
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)