

Curcumin (high purity)

Inhibitor of Lipoygenase and COX

Dual inhibitor of 5-lipoxygenase ($IC_{50}=8\mu M$) and cyclooxygenase (COX) ($IC_{50}=52\mu M$). Also inhibits glutathione S-transferase, induction of nitric oxide (NO) in activated macrophages and inhibits EGF-induced tyrosine phosphorylation of EGF receptors. Antioxidant. Inhibitor of NAD(P)H:quinone oxidoreductase 1 (NQO1), disrupting the binding of NQO1 to wild type p53 inducing ubiquitin-independent degradation of p53 and inhibits p53-mediated apoptosis in normal thymocytes and myeloid leukemic cells. Has been shown to inhibit NF- κ B, possibly through inhibition of JNK ($IC_{50}=10\mu M$), the COP9 signalosome kinase ($IC_{50}=10\mu M$) and p300/CBP.

Citations: 39

[View Online »](#)

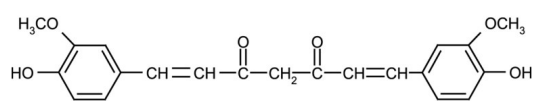
Ordering Information

[Order Online »](#)

ALX-350-028-M010	10mg
ALX-350-028-M050	50mg
ALX-350-028-M250	250mg

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	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1,7-bis(4-Hydroxy-3-methoxyphenyl)-1,6-heptadiene-3,5-dione, Diferuloylmethane
Appearance	Orange-yellow crystalline powder.
CAS	458-37-7
Couple Target	COX, GST, Lipoxygenase, NF-kappaB, NQO
Couple Type	Inhibitor
Formula	$C_{21}H_{20}O_6$
MI	14: 2673
MW	368.4
Purity	≥98% (HPLC)
Solubility	Soluble in acetic acid, 100% ethanol (at about 1mg/ml) or DMSO (25mg/ml).
Source	Isolated from turmeric (<i>Curcuma longa</i>).



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-
be@enzolifesciences.com](mailto:info-be@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)