

Xanthohumol

Prenylated flavonoid found in *Humulus lupulus*.

Ordering Information

[Order Online »](#)

LKT-X0254-M005	5mg
LKT-X0254-M010	10mg
LKT-X0254-M025	25mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

CAS 6754-58-1

Formula $C_{21}H_{22}O_5$

MW 354.4

Purity ≥98%

Solubility The solubility of xanthohumol in ethanol and DMF is approximately 3 mg/ml and approximately 2.5 mg/ml in DMSO. Xanthohumol is sparingly soluble in aqueous buffers.



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-be@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