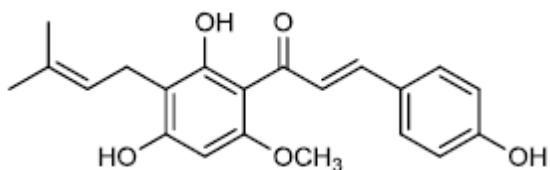


Xanthohumol

Cat. # AC-1008

Background: Binds to the N-terminal domain of valosin-containing protein (VCP), an essential protein for autophagosome maturation. Xanthohumol inhibits the function of VCP thereby impairing autophagosome maturation resulting in accumulation of the microtubule-associated protein 1 light chain 3-II (LC3-II)¹. Impairs prostate cancer cell growth and proliferation². Suppresses growth of pancreatic cancer cells via down-regulation of Notch1 signaling³. Cell permeable. Xanthohumol, a prenylated chalcone from hop, inhibits COX-1 and COX-2 activity and shows chemopreventive effects.

Application: Impairs autophagosome maturation



Product Information

Purity:	>98% by TLC
MW:	354.4
Formula:	C ₂₁ H ₂₂ O ₅
CAS No.	6754-58-1
Physical State:	Lyophilized yellow powder
Amount:	5mg, 25mg
Solubility:	35 mg/mL in DMSO or ethanol
Storage:	Store desiccated as supplied at -20°C for up to 3 years. Store solutions at -20°C for up to 1 month.

References

1. Y Sasazawa *et al.* *ACS Chem. Biol.* 2012 7:892
2. R Vene *et al.* *Mol. Med.* 2012 18:1292
3. S Kunnimalaiyaan *et al.* *Mol. Cancer Ther.* 2015 14:1395

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2013 LifeSensors, Inc. All Rights Reserved