

Xanthohumol

Catalog Number P021-10MG

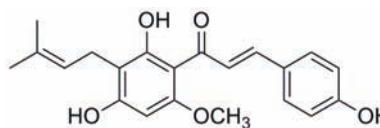
Catalog Number P021-50MG



ARBOR
ASSAYS

FEATURES

- Impairs prostate cancer cell growth
- Inhibits the function of VCP
- Cell Permeable



INTRODUCTION

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 and resulting in accumulation of the microtubule-associated protein 1 light chain 3-II (LC3-II)¹. Impairs prostate cancer cell growth and proliferation². Cell permeable.

| | |
|--------------------------|---|
| FORM: | Yellow solid |
| MOLECULAR WEIGHT: | 354.4 |
| STORAGE: | -20°C desiccated |
| FORMULA: | C ₂₁ H ₂₂ O ₅ |
| CAS NUMBER: | 6754-58-1 |
| OTHER NAMES: | (2E)-1-[2,4-Dihydroxy-6-methoxy-3-(3-methyl-2-buten-1-yl)phenyl]-3-(4-hydroxyphenyl)-2-propen-1-one |
| USES: | Soluble to 35 mg/mL in DMSO or ethanol |

REFERENCES:

1. Y Sasazawa et al. ACS Chem. Biol. 2012 7:892
2. R Vene et al. Mol. Med. 2012 18:1292

FOR RESEARCH USE ONLY

info@ArborAssays.com • www.ArborAssays.com

15Apr13