

## Curcumin

**Catalog #:** 27104

**Lot #:** 111121

**Size:** 1 g

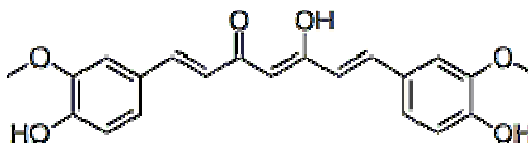
**Structure:**

**CAS Registry #:** 458-37-7

**Purity:** >99%

**Chemical Formula:** C<sub>21</sub>H<sub>20</sub>O<sub>6</sub>

**Molecular Weight:** 368.38



**Description:** Curcumin, also known as diferuloylmethane, natural yellow 3, or C.I. 75300, is the principal curcuminoid of the popular Indian spice turmeric, which is a small molecule, selective HAT inhibitor. It has shown to promote degradation of p300 and inhibits the acetyltransferases activity of purified p300 using either histone H3 or p53 as a substrate.

**Appearance:** Yellow powder.

**Solubility:** Soluble in acetone, ethanol and DMSO.

**Biological Activity:** Acetylation of histones H3 and H4 by p300/CBP was strongly inhibited by curcumin with an IC<sub>50</sub> 25 μM.

**Storage/Stability:** Store at or below -20 °C.

**Quality Control:** The purity was determined by HPLC analysis.

### Reference(s):

1. Marcu, MG. *et al.*, *Med. Chem.* 2006 Mar; **2(2)**: 169-174.
2. Kundu, TK. *et al.*, *Journal of Biological Chemistry* (2004); **279(49)**:51163-51171.