

---

## Curcumin

(Catalog No. M35000)

---

### Alternative Names

(E,E)-1,7-bis(4-Hydroxy-3-methoxyphenyl)-1,6-heptadiene-3,5-dione Diferuloylmethane  
Diferulylmethane

### Description

Curcumin is a yellow-orange dye obtained from tumeric, the powdered root of *Curcuma longa*; it is employed in the preparation of curcuma paper, the detection of boron, and possesses a vast number of biological targets. Curcumin has been shown to inhibit NF- $\kappa$ B (possibly thru inhibition of JNK; IC<sub>50</sub>=10  $\mu$ M), blocks amyloid peptide induced expression of: TNF- $\alpha$ , IL-1 $\beta$ , MCP-1, IL-8, and CCR5, regulates NOS2, inhibits IgE and Ag-induced degranulation of mast cells, inhibits 5-LO (5-lipoxygenase; IC<sub>50</sub>=8  $\mu$ M), Cox-2 (cyclooxygenase-2; IC<sub>50</sub>=52  $\mu$ M), the COP9 signalosome kinase (IC<sub>50</sub>=10  $\mu$ M), and p300/CBP. This wide range of activities makes Curcumin an extremely useful tool for developing novel analogous molecules and discovering new uses for this naturally occurring compound.

### CAS Number

458-37-7

### Chemical Formula

C<sub>21</sub>H<sub>20</sub>O<sub>6</sub>

### Molecular Weight

368.38

### Size

1 g

### Purity

≥95%

### Solubility

Soluble in water, ethanol (10 mg/ml), DMSO (25 mg/ml) and acetic acid

### Storage Temp

-20°C

### Ordering Information:

Products	Size	Cat. No.
Curcumin	1 g	M35000 -1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*