

Ellagic Acid (TBBD)

Catalog #: 27219

Lot #: 120608

Size: 100 mg

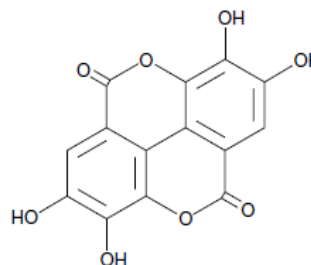
Structure:

CAS Registry #: 476-66-4

Purity: ≥95%

Chemical Formula: C₁₄H₆O₈

Molecular Weight: 302.2



Description: Ellagic acid, also known as TBBD, Gallogen, and Lagistase, blocks the methylation of arginine 17 of histone 3 (H3R17) by coactivator-associated arginine methyltransferase 1 (CARM1/PRMT4), without significantly altering histone acetylase or DNA methyltransferase activity. Ellagic acid can also alter cytochrome P450 activity, improve metabolism and clearance of xenobiotics, and alter immune function.

Appearance: A crystalline solid

Solubility: Soluble in DMSO, purged with an inert gas. Solubility in DMSO is ~0.14 mg/ml. Do not store aqueous solutions for more than one day.

Biological Activity: Ellagic acid was found to partially inhibit CARM1, even at an inhibitor concentration 10-fold above the IC₅₀ of 25 μM.

Storage/Stability: Store at or below -20°C for up to two years.

Quality Control: The purity was determined by HPLC analysis.

References:

1. Allen, C.T., *et al.* *Immunopharmacol. Immunotoxicol.* 2003; **25(3)**:409-422.
2. Herber, D., *et al.*, *Cancer Lett.* 2008; **269**:262-268.
3. Girish, C., *et al.*, *Fundam. Clin. Pharmacol.* 2008; **22**:623-632.
4. Selvi, B.R., *et al.*, *J. Biol. Chem.* 2010 Mar 5; **285(10)**: 7143-7152.
5. Paluszczak, J., *et al.*, *Toxicol. Lett.* 2010; **192(2)**:119-125.
6. Lea, M.A., *et al.*, *Anticancer Res.* 2010; **30(2)**:311-318.