

## Quercetin

**Catalog #:** 27214

**Lot #:** 120605

**Size:** 5 g

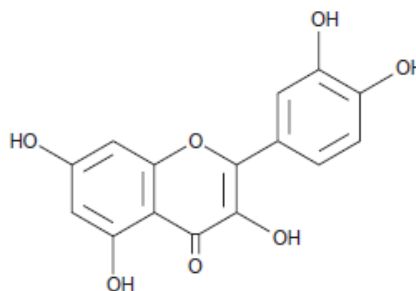
**Structure:**

**CAS Registry #:** 117-39-5

**Purity:** >95%

**Chemical Formula:** C<sub>15</sub>H<sub>10</sub>O<sub>7</sub>

**Molecular Weight:** 302.2



**Description:** Quercetin is an inhibitor of PDEs of both cAMP and cGMP. It is a flavonoid found in plant and fruit bark or rinds. It is estimated that normal dietary intake of Quercetin for humans is 0.1-0.2 mg/kg. Quercetin can induce renal adenomas in male rats when fed at 2,000 mg/kg.

**Appearance:** A crystalline solid

**Solubility:** Soluble in ethanol, DMSO and DMF, purged with an inert gas. Solubility in ethanol is 2 mg/ml and solubility in DMSO and DMF is 30 mg/ml. For maximum solubility in aqueous buffers, dissolve in DMSO and then dilute with the aqueous buffer of choice. Do not store aqueous solutions for more than one day.

**Biological Activity:** Quercetin has been shown to inhibit PDE 5A with an IC<sub>50</sub> value of 1.9 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. Gschwendt, M., *et al.*, *Biochem. Biophys. Res. Commun.* 1983; **117(2)**: 444-447.
2. Young, J.F., *et al.*, *Am. J. Clin. Nutr.* 1999; **69(1)**: 87-94.
3. Toxicology and carcinogenesis studies of quercetin in F344/N rats. National Toxicology Program, Technical Report, TR-409, 1-171. Retrieved August 4, 2004, from <http://ntp-server.niehs.nih.gov/htdocs/LT-studies/tr409.html>
4. Ruckstuhl, M., *et al.*, *Biochem. Pharmacol.* 1979; **28(4)**: 535-538.
5. Lang, D.R., *et al.*, *Biochim. Biophys. Acta.* 1974; **333(2)**: 180-186.
6. Lines, T.C., *et al.*, *Phytomedicine.* 2006 Mar; **13(4)**: 236-9.