
Epigallocatechin Gallate

(Catalog No. M21000)

Alternative Names

EGCG; (2R,3R)-2-(3,4,5-Trihydroxyphenyl)-3,4-dihydro-1[2H]-benzopyran-3,5,7-triol 3-(3,4,5-trihydroxybenzoate)

Description

A polyphenol flavonoid that displays antitumor and antioxidant properties. Inhibits telomerase and DNA methyltransferase (DNMT) and blocks the activation of EGF receptors and HER-2 receptors. Also potently and specifically inhibits chymotrypsin-like activity of the proteasome in vitro (IC₅₀ = 86-194 nM) and in vivo (IC₅₀ = 1 -10 μM).

CAS Number

989-51-5

Chemical Formula

C₂₂H₁₈O₁₁

Molecular Weight

458.37

Size

50 mg

Purity

≥98% by HPLC

Solubility

DMSO or EtOH

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
Epigallocatechin Gallate	50 mg	M21000-1

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