

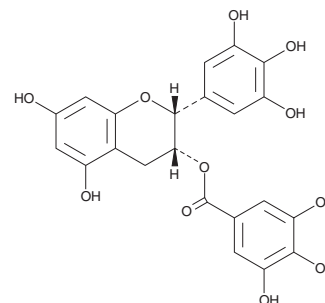
PRODUCT INFORMATION



(-)-Epigallocatechin Gallate

Item No. 70935

CAS Registry No.: 989-51-5
Formal Name: 3,4-dihydro-5,7-dihydroxy-2R-(3,4,5-trihydroxyphenyl)-2H-1-benzopyran-3R-yl-3,4,5-trihydroxy-benzoate
Synonym: EGCG
MF: C₂₂H₁₈O₁₁
FW: 458.4
Purity: ≥96%
UV/Vis.: λ_{max}: 276 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years
Item Origin: Plant/*Folium camelliae*



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

(-)-Epigallocatechin gallate (EGCG) is supplied as a crystalline solid. A stock solution may be made by dissolving the EGCG in an organic solvent purged with an inert gas. EGCG is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of EGCG in these solvents is approximately 20, 25, and 30 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of EGCG can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of EGCG in PBS (pH 7.2) is approximately 25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

EGCG is a phenol that has been found in a variety of plants including green and black tea plants and has diverse biological activities.¹⁻⁶ It is lytic against *T. cruzi* amastigotes, with a 50% bactericidal concentration (MBC₅₀) value of 0.53 pM.¹ EGCG decreases oxidation of deoxyguanosine and tyrosine nitration (IC_{50s} = 0.25 and 0.11 mM, respectively) and inhibits macrophage-mediated LDL oxidation *in vitro*.^{2,3} EGCG increases LDL receptor binding activity and lowers cholesterol levels in HepG2 cells *via* increased cholesterol efflux.⁴ It inhibits heregulin-β1-induced activation of ErbB2 and ErbB3 as well as migration and invasion of MCF-7 breast cancer cells.⁵ *In vivo*, EGCG reduces food intake and body weight, serum levels of insulin, leptin, testosterone, and growth hormone, and the size of the ventral prostate, testis, liver, and spleen in male Sprague Dawley, lean Zucker, and obese Zucker rats when administered at doses of 82, 81, and 92 mg/kg, respectively.⁶

References

1. Paveto, C., Güida, M.C., Esteva, M.I., *et al.* *Antimicrob. Agents Chemother.* **48(1)**, 69-74 (2004).
2. Fiala, E.S., Sodum, R.S., Bhattacharya, M., *et al.* *Experientia* **52(9)**, 922-926 (1996)
3. Yoshida, H., Ishikawa, T., Hosoai, H., *et al.* *Biochem. Pharmacol.* **58(11)**, 1695-1703 (1999)
4. Bursill, C.A. and Roach, P.D. *J. Agric. Food Chem.* **54(5)**, 1621-1626 (2006).
5. Kushima, Y., Iida, K., Nagaoka, Y., *et al.* *Biol. Pharm. Bull.* **32(5)**, 899-904 (2009).
6. Kao, Y.H., Hiipakka, R.A., and Liao, S. *Endocrinol.* **141(3)**, 980-987 (2000).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/15/2019

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM