# **PRODUCT** INFORMATION



## **Echinacoside**

Item No. 29927

CAS Registry No.:	82854-37-3	
Formal Name:	2-(3,4-dihydroxyphenyl)ethyl	ОН
	O-6-deoxy-α-L-mannopyranosyl-	
	(1→3)-O-[β-D-glucopyranosyl-	HONOH
	$(1 \rightarrow 6)$ ]-β-D-glucopyranoside,	
	4-[3-(3,4-dihydroxyphenyl)-2-propenoate]	
Synonyms:	Kusaginin, NSC 603831, trans-Acteoside,	
	trans-Verbascoside	
MF:	$C_{35}H_{46}O_{20}$	
FW:	786.7	ОН ОН
Purity:	≥95%	
UV/Vis.:	λ <sub>max</sub> : 219, 336 nm	HOʻ
Supplied as:	A crystalline solid	НО
Storage:	-20°C	ОН
Stability:	≥4 years	
Item Origin:	Plant/Cistanches herba	

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

#### Laboratory Procedures

Echinacoside is supplied as a crystalline solid. A stock solution may be made by dissolving the echinacoside in the solvent of choice, which should be purged with an inert gas. Echinacoside is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of echinacoside in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Echinacoside is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, echinacoside should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Echinacoside has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

#### Description

Echinacoside is a phenylethanoid glycoside that has been found in Echinacea and has diverse biological activities.<sup>1-4</sup> It scavenges 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) radicals  $(EC_{50} = 6.6 \ \mu M$  in a cell-free assay).<sup>1</sup> Echinacoside (10  $\mu g/ml$ ) inhibits hydrogen peroxide-induced production of reactive oxygen species (ROS) and apoptosis in PC12 cells.<sup>2</sup> It reduces increases in plasma alanine aminotransferase (ALT), TNF- $\alpha$ , IL-1 $\beta$ , IL-16, and IL-10 levels and hepatocyte apoptosis induced by LPS/D-galactosamine in a mouse model of acute liver injury when administered at a dose of 60 mg/kg.<sup>3</sup> Echinacoside (20 mg/kg per day for 15 days) prevents loss of dopaminergic neurons in the substantia nigra pars compacta and decreases in striatal dopamine and 3,4-dihydroxyphenylacetic acid (DOPAC) levels in a mouse model of MPTP-induced Parkinson's disease.<sup>4</sup> It also improves motor performance in the rotarod test in the same model.

### References

- 1. Pellati, F., Benvenuti, S., Magro, L., et al. J. Pharm. Biomed. Anal. 35(2), 289-301 (2004).
- 2. Kuang, R., Sun, Y., Yuan, W., et al. Planta Med. 75(14), 1499-1504 (2009).
- 3. Li, X., Gou, C., Yang, H., et al. Scand. J. Gastroenterol. 49(8), 993-1000 (2014).
- 4. Geng, X., Tian, X., Tu, P., et al. Eur. J. Pharmacol. 564(1-3), 66-74 (2007).

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

#### SAFFTY 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.

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, 10/31/2022

## 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