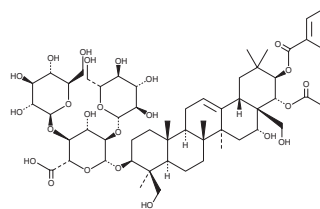


PRODUCT INFORMATION

Aescin

Item No. 28815

Synonym:	Escin
MF:	$C_{55}H_{86}O_{24}$ (for β -Aescin)
FW:	1,131.3
Purity:	$\geq 95\%$
Supplied as:	A solid
Storage:	-20°C
Stability:	≥ 2 years
Item Origin:	Plant/Horse chestnut



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

Laboratory Procedures

Aescin is supplied as a solid. A stock solution may be made by dissolving the aescin in the solvent of choice, which should be purged with an inert gas. Aescin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of aescin in these solvents is approximately 25 and 20 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 aescin can be prepared by directly dissolving the solid in aqueous buffers. The solubility of aescin in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Aescin is a mixture of aescin congeners and isomers, which are saponins that have been found in *A. hippocastanum*.^{1,2}

References

1. Sirtori, C.R. Aescin: Pharmacology, pharmacokinetics and therapeutic profile. *Pharmacol. Res.* **44**(3), 183-193 (2001).
2. Savarino, P., Colson, E., André, J., *et al.* Horse chestnut saponins-escins, isoescins, transescins, and desacylescins. *Molecules* **28**(5), 2087 (2023).

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, 09/06/2023

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