

PRODUCT INFORMATION



Cholestyramine

Item No. 37827

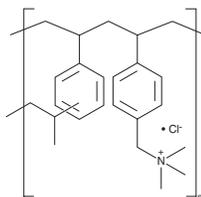
CAS Registry No.: 11041-12-6

Synonym: MK-135

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Description

Cholestyramine is a bile acid sequestrant.¹ It binds to bile acids in the intestine and inhibits their absorption, which increases the synthesis of bile acids from cholesterol in the liver, thereby lowering serum LDL levels. Cholestyramine reduces serum cholesterol levels and increases bile acid excretion in the feces without changing the bile acid pool size in rabbits.² Formulations containing cholestyramine have been used as adjuncts in the treatment of primary hypercholesterolemia.

References

1. Scaldaferrri, F., Pizzoferrato, M., Ponziani, F.R., *et al.* Use and indications of cholestyramine and bile acid sequestrants. *Intern. Emerg. Med.* **8(3)**, 205-210 (2013).
2. Qiu, D.N., Shang, Q., Sun, D.Y., *et al.* Cholestyramine alters bile acid amounts and the expression of cholesterol-related genes in rabbit intestinal and hepatic tissues. *J. Dig. Dis.* **18(2)**, 107-114 (2017).

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, 05/30/2024

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