

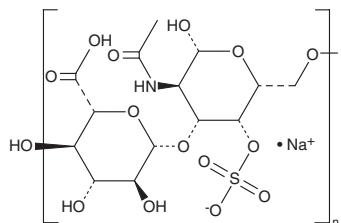
PRODUCT INFORMATION



Chondroitin Sulfate (sodium salt)

Item No. 37801

CAS Registry No.: 9082-07-9
Formal Name: chondroitin, hydrogen sulfate, monosodium salt
Synonyms: Chondroitin-4-sulfate, CSA
MF: $(C_{14}H_{22}NO_{15}S)_n \cdot Na$
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.

Laboratory Procedures

Chondroitin sulfate (sodium salt) is supplied as a solid. Aqueous solutions of chondroitin sulfate (sodium salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of chondroitin sulfate (sodium salt) in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Chondroitin sulfate is a glycosaminoglycan (GAG).^{1,2} It contains a GAG-protein linker region and repeating disaccharide subunits composed of glucuronic acid (GlcA) and N-acetylgalactosamine (GalNAc) that are often sulfonated. Chondroitin sulfate chains of various lengths covalently attach to the surface of proteins via a serine residue in the GAG-protein linker region, forming chondroitin sulfate proteoglycans (CSPGs).² CSPGs participate in the regulation of many biological activities, such as cell adhesion and migration, organization of brain extracellular matrix (ECM), and the anti-inflammatory response.²⁻⁴

References

1. Kwok, J.C.F., Warren, P., and Fawcett, J.W. Chondroitin sulfate: A key molecule in the brain matrix. *Int. J. Biochem. Cell Biol.* **44**(4), 582-586 (2012).
2. Weisner, J., Landel, I., Reintjes, C., *et al.* Preclinical efficacy of covalent-allosteric AKT inhibitor borussertib in combination with trametinib in KRAS-mutant pancreatic and colorectal cancer. *Cancer Res.* **79**(9), 2367-2378 (2019).
3. Mikami, T. and Kitagawa, H. Biosynthesis and function of chondroitin sulfate. *Biochim. Biophys. Acta* **1830**(10), 4719-4733 (2013).
4. Yamaguchi, Y. Lecticans: Organizers of the brain extracellular matrix. *Cell. Mol. Life Sci.* **57**(2), 276-289 (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, 05/08/2026

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