

Certificate of Analysis



Product Name: Sodium Hyaluronate, 2000 kDa

Product Number: H-2000

Lot Number: PHI4004B

Product Appearance: White or almost white, very hygroscopic fibrous aggregate

Average MW: 2,200 kDa by Size Exclusion Chromatography

Polydispersity Index: 1.3

IR Spectrum: Complies to the Ph. Eur. Reference spectrum of sodium hyaluronate.

Intrinsic Viscosity: 2.40 m³/kg

Proteins: Not detected, <0.01%

Chloride: <0.124%

Iron: <5 ppm

Loss on Drying: 12.6%

Bacterial Endotoxins: <0.020 I.U./mg

Sodium Hyaluronate content / dried product: 98.0%

Passes All Functional Specifications

A handwritten signature in black ink, appearing to be 'Colin Ferguson', written over a horizontal line.

4/16/2021

Colin Ferguson
Quality Manager

**Sodium Hyaluronate is manufactured by HTL and distributed by Echelon.
For research use only. Not intended for diagnostic or therapeutic use.**

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. This product is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development. or health or safety of any uses of the product.