



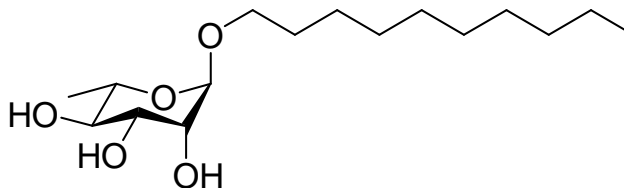
Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only
Not intended or approved for
diagnostic or therapeutic use.

Product Name: Decyl Rhamnoside (decyl- α -L-rhamnopyranoside) >95%

Product Number: RL10-1



Molecular Formula: C₁₆H₃₂O₅ **MW:** 304.42 **CAS:** N/A

Solubility: Soluble in water ~18mg/L, soluble in most organic solvents.

Storage: room temperature or below

Critical Micelle Concentration (CMC): ~344 μ M

Surface Tension (γ CMC): ~29 mN/m

pH = 6 – 8 in water

Estimated HLB: 11.8 (classified as a water-soluble emulsifier and wetting agent)

Field of Interest: Decyl rhamnoside surfactants are suitable for use in cosmetics, including in anti-aging creams, soaps, and other skin or facial treatments. Furthermore, these materials may have anti-bacterial properties, anti-zoosporic properties, and anti-fungal. Agriculture uses are known, as well as bioremediation, and enhanced oil recovery projects.

General: Decyl rhamnoside (like many of its natural rhamnose-based derivatives) is a good foaming and wetting agent, with the ability to increase the aqueous solubility of hydrophobic compounds, making them excellent solubilizing and emulsifying agents for diverse applications.

Produced in collaboration with GlycoSurf, LLC. (www.glycosurf.com)

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. Decyl Rhamnoside 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.

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at busdev@echelon-inc.com.

TDS RL10-1 Rev: 2 (04/07/2015)