



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	Cat#	MW
HA850-TexasRed	H-700R	680 kDa
HA300-TexasRed	H-250R	240 kDa
HA30-TexasRed	H-025R	24 kDa

**Solubility:** H<sub>2</sub>O

**Storage:** - 20 °C

**Field of Interest:** Cell Biology, Biomaterials, Tissue Engineering, Drug Delivery

**Absorption/Emission( $\lambda_{\max}/\text{nm}$ ):** 589/615

#### References:

1. Hyaluronic acid based scaffolds for tissue engineering --- A review, Carbohydrate Research 2013, 92(2):1262-1279
2. Controlled chemical modification of hyaluronic acid: synthesis, applications, and biodegradation of hydrazide derivatives, Journal of Controlled Release, 1998, 53(1-3): 93-103
3. Self-assembled hyaluronic acid nanoparticles for active tumor targeting, Biomaterials, 2010, 31, 106–114

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