



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: Oral Formulation Vehicle

Product Number: B-0706

Storage: 4° C

Suggested Use: This buffer is intended to create aqueous suspensions of organic soluble compounds for oral administration in nonhuman research subjects. The vehicle is formulated with a cellulose compound not metabolized by laboratory rodents.