



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: SphingoBeads™
Catalog Number: S-6100

QUANTITIES AVAILABLE: 1 ml/vial

SPECIFICATIONS: One ml of agarose beads containing ~10 nanomoles of bound sphingosine. Beads are in PBS buffer. Each order of Sphingo-Beads™ includes a small amount of control beads.

**STORAGE
RECOMMENDATIONS:** Store at 2-8°C. Do not freeze.

SUGGESTED USAGE: lipid-protein pull-down.