



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.

## SHIP2 Reaction Buffer

**Product Number:** K-S2RB

**Storage:** Store at 4°C

**Description:** SHIP2 Enzyme Reaction Buffer. Use for measuring inositol polyphosphate 5-phosphatase activity. Echelon's SHIP2 reaction buffer recipe is proprietary.

**Size:** 10 mL

**Suggested Use:** Use this SHIP2 reaction buffer for measuring inositol polyphosphate 5-phosphatase activity or for SHIP2 inhibitor study. This buffer is ready for use. Dilute SHIP2 enzyme and substrate directly in this buffer. This buffer also works for Schizosaccharomyces Pombe Synaptojanin.

### Suggested Reaction Conditions:

- Malachite Green Assay (product # K-1500)  
25 µL reaction with 120 µM PI(3,4,5)P<sub>3</sub> substrate (product # P-3908, 3 nmol per reaction) and 5-20 nM SHIP2 (product # E-1000) in SHIP2 reaction buffer for 30 – 60 min at 37°C.
- 5'-PtdIns(3,4,5)P<sub>3</sub> Phosphatase Activity Fluorescence Polarization Assay (product # K-1400)  
2 µM PI(3,4,5)P<sub>3</sub> substrate (part # K-1403L) and 5-10 nM SHIP2 (product # E-1000) in SHIP2 reaction buffer for 30 – 60 min at 37°C.

**Notes:** Optimization of the phosphatase reactions by enzyme and PI(3,4,5)P<sub>3</sub> substrate titration, and reaction time may be required.