



P.O. Box 58537  
Salt Lake City, Ut 84108  
Telephone 801-588-0455  
Fax 801-588-0497  
Email: echelon@echelon-inc.com  
http://www.echelon-inc.com

## Certificate of Analysis

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

**Product Name:** Spingosine 1 Phosphate Fluorescein

**Product Number:** S-200F

**Lot Number:** E00138-78

**Product Description:** Yellow powder

**Quality Assay:**  $^1\text{H}$  and  $^{31}\text{P}$  NMR, Thin Layer Chromatography

### Quality Parameters

- A. NMR, Spectra matches control spectra
- B. Thin Layer Chromatography

### Passes

✓  
✓

Passes All Functional Specifications



Date 01/20/2010

Lee Crosby  
Quality Manager