

Catalog #: 81011

**AnaSpec, Inc.**

## Analytical Data Sheet

---

**Product Name:** 6-JOE,SE(6-carboxy-4,5-dichloro-2,7-dimethoxyfluorescein,succinimidyl ester

**Lot No.:** CM6094

**Weight:** 5mg

**Appearance:** Red solid

**Formula:**

**Molecular Weight:**

**Melting Point:**

**Loss on Drying:**

**Free Amino Acids:**

**HPLC**

**Additional Instructions:** For best results, pre-swell resin with DCM for 10 minutes.

---

**Remarks: Not for Human Use. Research Purposes Only.**

**Release By:** \_\_\_\_\_ **Date:** \_\_\_\_\_  
Quality Control