

Catalog #:AS-81007

# AnaSpec, Inc.

## Analytical Data Sheet

---

<b>Product Name:</b>	<b>5-FAM,SE (5-carboxyfluorescein,succinimidyl ester)</b>
<b>Lot No.:</b>	CM1653
<b>Weight:</b>	10mg
<b>Formula:</b>	$C_{25}H_{15}NO_9$
<b>Molecular Weight:</b>	473.39
<b>Appearance:</b>	Orange solid
<b>Purity:</b>	
<b>TLC:</b>	>90%
<b>HPLC:</b>	>90% purity

---

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

**Release By:** \_\_\_\_\_ **Date:** \_\_\_\_\_

*Quality Control*