

Catalog #: 81004

AnaSpec, Inc.

Analytical Data Sheet

Product Name: 6-FAM (6-carboxyfluorescein)

Lot No.: CM2096

Weight: 100 mg

Formula: $C_{21}H_{12}O_7$

Molecular Weight: 376.32

Appearance: Red solid

Purity:

TLC: >95 % purity

Remarks: Not for Human Use. Research Purposes Only.

Release By: _____ **Date:** _____

Quality Control