

Catalog #:AS-60438-F1

# AnaSpec, Inc.

## Analytical Data Sheet

---

**Product Name:** Fmoc-trans-4-Fluoro-Pro-OH

**Lot No.:** AA34274

**Weight:** 1g

**Formula:** C<sub>20</sub>H<sub>18</sub>FNO<sub>4</sub>

**Molecular Weight:** 355.4

**Appearance:** Off-white solid

**HPLC:** 98% purity

**Solubility:** 1mmol in 1ml DMF

**Storage:** Keep in a cool & dry place

---

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

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

*Quality Control*