

---

## CERTIFICATE OF ANALYSIS

**Catalog Number** C020  
**Product Name** Pluronic® F-127  
**CAS Number** 9003-11-6  
**Molecular Weight** ~12500  
**Appearance** white solid  
**Lot Number**

	LOT DATA	SPECIFICATION
<b>Absorption<sup>1</sup></b> Optical Density at 280 nm	0.013	<0.10

1. Measure at 10 mg/mL in E-pure water.