



## Product Data Sheet

---

Product Name:	Ninhydrin Test Kit for Resin		
Catalog Number:	AS-25241	Lot Number:	See label on vial
Description:	This kit consists of three bottles (10 ml each) of reagents. It is designed to check if the coupling of a protected amino acid to the amino group of the resin is complete.		
Storage:	2 - 8 °C		
Test Protocol:	<ol style="list-style-type: none"><li>1. Place 20-30 beads of resin in a 12x75 test tube.</li><li>2. Add 2 drops of reagent 1, 2 and 3 to the test tube.</li><li>3. Heat the test tube in a beating block at approximately 120 °C for 3 minutes.</li><li>4. Data Analysis<ol style="list-style-type: none"><li>4.1 Yellowish color: Complete coupling</li><li>4.2 Blue color: Incomplete coupling</li></ol></li></ol>		

*For Research Use Only*