



# Recombinant Swine IL-10 ELISA Standard

## PRODUCT ANALYSIS SHEET

---

<b>Catalog Number:</b>	SD064
<b>Lot Number:</b>	See product label
<b>Quantity:</b>	Refer to vial label for mass and reconstitution volume.
<b>Formulation:</b>	Lyophilized in a solution containing PBS, 3% BSA, stabilizers and 15 mM sodium azide (Caution: sodium azide is a poisonous and hazardous substance. Handle with care and dispose of properly.)
<b>Storage:</b>	The lyophilized preparation should be stored at 2 to 8°C. Once reconstituted, store the product in working aliquots at $\leq -80^{\circ}\text{C}$ . This product may be subjected to a single freeze/thaw without significant loss of reactivity in ELISA.
<b>Expiration Date:</b>	Lyophilized material expires one year from date of receipt when stored as instructed.

**This product is for research use only. Not for use in diagnostic procedures.**

[www.invitrogen.com](http://www.invitrogen.com)

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PISD064 (Rev 01/09)

(Page 1 of 1) DCC-09-0057

**Important Licensing Information** - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.