

# invitrogen™ Human Tumor Necrosis Factor Alpha ( TNF- $\alpha$ ) Control Set

---

<b>Catalog Number:</b>	CS003
<b>Lot Number:</b>	See product label
<b>Quantity:</b>	1 vial each Control 1 and Control 2. See vial label for concentrations.
<b>Formulation:</b>	Lyophilized.
<b>Reconstitution Recommendations:</b>	Reconstitute each vial with 2.0 mL distilled water or other appropriate matrix. Allow the controls to rehydrate for 10 minutes, then swirl gently to mix.
<b>Storage:</b>	Store lyophilized human TNF- $\alpha$ Controls at 2-8°C, preferably desiccated. Upon reconstitution, aliquot into polypropylene tubes and store at -80°C. Do not subject to more than one freeze-thaw cycle.
<b>Expiration Date:</b>	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 - Carlsbad - CA 92008 - Tel: 800.955.6288 - E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)