



## Human Heart Mitochondria WB Control

<b>CATALOG #:</b>	438505
<b>COMPONENTS:</b>	50 µg human heart mitochondria
<b>APPLICATIONS:</b>	To be used as a positive control for Western blotting
<b>CONCENTRATION:</b>	1 µg/µL in SDS-PAGE sample buffer (2% SDS, 60 mM Tris, pH 6.8, with HCl, 5% glycerol, 0.02 mg/mL Bromophenol Blue), 50 mM DTT.
<b>SUGGESTED WORKING CONCENTRATION:</b>	0.5 – 5.0 µg/lane for Western blotting
<b>STORAGE CONDITIONS:</b>	Shipped at 4°C. Store at -20°C after receipt or -80°C for long term.
<b>COUNTRY OF ORIGIN:</b>	USA
<b>ADDITIONAL INFO:</b>	<u>Human Tissue Handling Risks and Precautions:</u> These samples were prepared post-mortem from a donor whose serology report was negative for Hepatitis C antibody, HIV-1 & 2, Hepatitis B Surface antigen. However, as with all human derived material, standard lab practices to prevent direct contact should be used.

---

### For Research Use Only



Manufactured exclusively for Invitrogen by MitoSciences, Inc.

[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)

PI438505

(Rev 10/08) DCC-08-1089

**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.