

## Certificate of Analysis

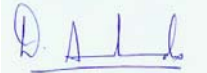
Product: Mouse Laminin-I  
PathClear®  
Catalog Number: 3400-010-02  
Size: 1 ml (1 mg/ml)  
Storage: Store at -80°C

Description: **SPECIFICATION CERTIFICATE**

Test/Method	Specification	Results
Sterility Testing (USP XXIII)	Sterile for Bacteria and Fungi	PASS
Endotoxin	≤20 EU/ml by LAL Assay	PASS
Purity	> 90% by SDS-PAGE	PASS
Functional Assays	Promotes attachment and formation of neuronal filaments in NG108-15 cells.	PASS
MAP Tests	<i>Mycoplasma pulmonis</i> <i>E. cuniculi</i> Viruses Tested: MHV, TMEV, MMV, PVM, MVM HANTAAN, SENDAI, MTV, MCMV ECTROMELIA, LCMV, LDEV POLYOMA, K Virus, REO(1,3) MPV, EDIM, MAV(1,2)	PASS PASS PASS
PCR Tests	<i>Mycoplasma</i> spp. Viruses Tested: LDEV, SENDAI, MHV, PVM, MVM MPV(1,2,3), NOROVIRUS, REO3 EDIM, ECTROMELIA, LCMV, K, MTV POLYOMA, HANTAAN, MAD(1,2) MCMV, KILHAM's, TOOLAN's, PARVO RCMV, CORONA, RMV, SEOUL SIALODACRYOADENITIS, TMEV, TMELV	PASS PASS

Quality Control:

Release: Lot Number: 16064F8  
Tested By: sfoster  
Retest Date: 6/27/2009  
Date: 06/30/2008

Approval: 

**Production Manager**

Trevigen, Inc.  
8405 Helgerman Ct.  
Gaithersburg, MD 20877  
301-216-2800 • 1-800-873-8443  
FAX: 301 560 4973  
e-mail: info@trevigen.com  
www.trevigen.com