



## RELATED PRODUCTS

Product	Catalog Code
LPS-EB Ultrapure	tlrl-pelps
pUNO-hTLR4A	puno-htrl4a
pUNO-mTLR4	puno-mtrl4
pDUO2-hMD2/CD14	pduo2-hmd2cd14
pDUO2-mMD2/CD14	pduo2-mmd2cd14
293/hTLR4A-MD2-CD14	293-htrl4amd2cd14
pNiFty-Luc (Amp <sup>R</sup> )	pnifty-luc
pNiFty-SEAP (Amp <sup>R</sup> )	pnifty-seap

---

### TECHNICAL SUPPORT

Toll free (US): 888-457-5873  
Outside US: (+1) 858-457-5873  
Europe: +33 562-71-69-39  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)  
Website: [www.invivogen.com](http://www.invivogen.com)



3950 Sorrento Valley Blvd. Suite A  
San Diego, CA 92121 - USA