



**Signosis, Inc.**

BioSignal Capture

528 Weddell Dr. Suite 5  
Sunnyvale, CA 94089, USA  
Tel: 408-747-0771  
Email: [info@signosisinc.com](mailto:info@signosisinc.com)  
Website: [www.signosisinc.com](http://www.signosisinc.com)

## New Account Form

**FAX BACK TO: 408-864-2182**

Institution / Company Name: \_\_\_\_\_

Purchasing Contact: \_\_\_\_\_ Tel: \_\_\_\_\_

**Bill To (Accts Payable) Information:**

Bill-To Contact: \_\_\_\_\_

Address: \_\_\_\_\_

\_\_\_\_\_

\_\_\_\_\_

Tel. \_\_\_\_\_

Fax: \_\_\_\_\_

Email: \_\_\_\_\_

**Ship To Information:**

Ship-To Contact: \_\_\_\_\_

Address: \_\_\_\_\_

\_\_\_\_\_

\_\_\_\_\_

Tel. \_\_\_\_\_

Fax: \_\_\_\_\_

Email: \_\_\_\_\_