





570-5600 Parkwood Way  
Richmond, BC V6V2M2  
CANADA  
Tel: 1-(604)-232-4600  
Fax: 1-(604)-232-4601  
orders@signalchem.com

Catalogue #	Description	Quantity	Unit Price	Total

**Total Order Amount:**

**Special Shipping Considerations / Additional Information:** ( if required )

**Promotion Code:**

For Internal Use Only	
CU:	TW:
BX:	DI:
FE:	
IV:	
DC:	