

ORDER FORM

FAX TO:

19520 Amaranth Drive
Germantown, Maryland 20874, USA

301-515-6983 (FAX)
301-515-6982 (Tel.)

PO No. _____

Promotion Code: _____

Shipping Address		Billing Address <input type="checkbox"/> Check Here if Same As Shipping Address	
Last Name		Attn.	
First Name		Street	
Department		Building/room	
Institution		City	
Street		State	
Building/room		Zip	
City		Country	
State		Tel	
Zip		Fax	
Country		Visa Card#	
E-mail		Master Card#	
Tel		Expiration Date	Month: Year:
Fax			

Cat. No.	Description	Price \$	Qty	Amount \$

_____ (Signature)

_____ (Date)