

CREDIT APPLICATION

Please submit your application to Orders@arborassays.com

Legal Name of Business:	Trade Name of Business (if different):	
Bill to Address:	Ship to Address:	
Phone #:	Fax #:	
Faraily	Condit Line De sweete de	
Email:	Credit Line Requested:	
Business Type:		
[] Partnership [] Limited Partr		
[] LLC [] Other (please	e specify)	
Federal ID#: Number of years in business:		
(If tax exempt, attach certificate)		
Accounts Payable Information:		
Name of AP Contact:	Title:	
Address:	Email:	
Phone #:	Fax #:	
Bank Information:		
Bank Name:	Account #:	
	Account Type:	
Bank Address:	Contact:	
	E-mail:	
Phone #:	Fax #:	
FIIOTIE #.	FdX #.	



CREDIT APPLICATION

Credit References:

Company: Contact: Phone #: Address:	Email: Fax #:
Company: Contact: Phone #: Address:	Email: Fax #:
Company: Contact: Phone #: Address:	Email: Fax #:
Terms: Payment in U.S. dollars (\$ I certify that the above information is Signature (Name and Title, please R	s correct and agree to the terms shown. (Date)