

SHIPPING CONDITIONS

VALID UNTIL DECEMBER 31st, 2017

- This price list supersedes all previous lists.
- Please mention your P.O. number, quotation number (for custom or bulk orders), as well as all applicable promotional codes on every order.
- Administrative costs will be charged for re-invoicing, if necessary, due to incorrect or incomplete information on the order form.
- Transport costs will be invoiced according to the table below or as priced in your quotation.
- Shipping fees are as follows for packages up to 10 lbs*:

		Standard Fee	Blue Ice Fee	Dry Ice Fee
US	Next-day delivery	\$35	\$10	\$20
	Two-day delivery	\$20		
Canada	Next-day delivery	\$60	\$10	\$20
International	Next-day delivery	\$60	\$20	\$80

*Additional fees may apply for hazardous chemicals

- All prices are exclusive of duties, taxes, customs and other fees.
- We are constantly improving our products and services. For this reason, all prices and specifications are subject to change without prior notice.

