

Important Note on Hazardous Chemical Shipping Charges

The U.S. Department of Transportation requires that 2-mercaptoethanol (Bio-Rad catalog number 161-0710) be packaged in specially approved boxes separate from other chemical goods. In addition, many carriers apply an extra charge for shipping hazardous chemicals. Because of this, special handling and shipping charges have been applied to your order. Consolidation of future orders will help minimize these charges in the future.