

**IBMX**

IBMX - Phosphodiesterase inhibitor - 40 µl

Product reference :

**62AMXADA**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 04A  
- Conditioning buffer \_\_\_\_\_ DMSO  
- Appearance \_\_\_\_\_ Liquid form  
- Storage conditions \_\_\_\_\_ 4 °C  
- Expiration date \_\_\_\_\_ 2023.03.19  
- Active Moiety Concentration (0) \_\_\_\_\_ 500 mM

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (0)	40.00 µl	222

*Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)*