

31 August 2011

Your Order Number:
Item:

Dear

Thank you for your recent order. As a manufacturer and supplier of research chemicals we are required to maintain strict controls over the supply of our chemicals and equipment. Your order contains an item which requires a declaration and/or licence. Below is a form which needs to be completed and faxed back to us before we can continue processing this product. Other items on your order not requiring a declaration will be despatched as per normal.

If your completed form is not returned to us within 10 working days the item will be cancelled.

Please fax back all forms to **01747 833569**

Please contact us if you have any questions regarding the forms as we want to process your order with the minimum delay. We can be contacted by telephone on 01747 833311

Sigma-Aldrich Customer Service

End Use Statement
TYPE OR PRINT CLEARLY

Our Ref: _____

Your PO Ref: _____

Product No: _____

Pack Quantity: _____

Product Description: _____

Company/Organisation Name: _____

Department Name: _____

Address: _____

(Post Office Box numbers are unacceptable)

Country: _____ **Postal Code:** _____

Telephone Number: _____ **Fax Number:** _____

End User: _____

(Full name please, no initials)

Detailed end usage. Please be specific, using common language where possible, and define any technical terms or jargon used. General statements such as “for research only” are not sufficient.

Are the goods going to be exported by yourself? *(Please tick)* YES NO

Are the goods going to be resold? *(Please tick)* YES NO

If the goods are to be resold, is the final customer exporting the goods? YES NO

I certify that, if none of the above then, I am the end user of the product(s) listed above and can therefore confirm that this product will NOT be: -

- a) Used for the development or manufacture of chemical or biological weapons
- b) Used for the development or manufacture of controlled drugs, as specified in “The Misuse of Drugs Act 1971” and its associated regulations
- c) Used directly, or in the formulation of preparations, for human or veterinary purposes.
- d) Used for any other purposes restricted by the conditions of sale shown in the current catalogue.

Signature: _____ **Date:** _____

Print Name: _____ **Job Title:** _____