

Export/Import of Controlled Substances Overseas



Why can it take so long to have my reference standard delivered overseas?

Delivery time of controlled reference materials requires the careful navigation of governmental agencies of both the country of origin and country of import. Several processes are in place to document the transfer of goods and additional steps are needed if the compound is recognized as an internationally or locally controlled substance.

Cayman has a dedicated regulatory specialist who processes and tracks these orders to help secure all necessary permits to facilitate timely delivery. Our international distributor network also helps provide local country support with customs clearance and delivery.

Timeline for the Export/Import of Controlled Substances



Export Procedures for US-Controlled Substances

- A foreign import certificate or 'Permit to Import' must be sent to Cayman's International US Drug Enforcement Agency (DEA) Order Processing Center.
- If the product is not considered a controlled substance in the importing country, then a letter of no objection issued by the appropriate competent national authority must be sent to Cayman.
- Once the import permit for an order is received, Cayman will file for an export permit from the DEA. This permit is typically issued within three to four weeks after the permit application is submitted.

International Narcotic Control Board (INCB) Resources

- Yellow and Green Lists of compounds under international control
- Status of estimates of narcotic drugs listed in Schedules I and II
- Assessment request forms for compounds that are not listed

Visit www.caymanchem.com/deliverytimes for links to these resources

IMPORT PERMIT CHECKLIST

- ☒ Determine order quantity is within your country's allotment
- ☒ Ensure the Permit to Import has at least 60 days before expiry
- ☒ Email permit to Cayman at international@caymanchem.com

Export Procedures for Items Not Controlled in the US

- The requisite customs import documents must be sent to Cayman so they can accompany the shipment.
- Export permits are not required for items not controlled in the US or for specific compounds in solution granted a US DEA exemption.

Shipping

- The compound is shipped only when both the import and export permits are granted.
- Delivery can take place in as little as two to three days depending on the customs arrangements of the importing country.

Learn more about the procedures and estimated timeline for the export/import of controlled substances by visiting www.caymanchem.com/deliverytimes

To find your local distributor, contact Keith Williams at kwilliams@caymanchem.com

If questions arise along the ordering process, contact Customer Service at custserv@caymanchem.com

Cayman Chemical Company
Attn: Int'l DEA Order Processing
1180 E Ellsworth Road
Ann Arbor, MI 48108
USA
Phone: (734) 975-3888

DEA Exempt Solutions

We make every attempt to offer exempt preparations as soon as possible. Some exempt preparations will have "exempt preparation" in the item name, while others will have "CRM" in the name. Not all CRM items are exempt preparations. You can use the search facets on the "Product & Attribute Search" on our Forensic page to filter your search by DEA Schedule.