



1360 Willow Road, Suite 205  
Menlo Park, CA 94025-1516  
Phone 650-322-1360

## MATERIAL SAFETY DATA SHEET

### PRODUCT IDENTIFICATION

Product Name:	Human IgG Calibrators; Q Assay
Catalogue Number:	18-1030
Unit Size:	Box with nine vials, each with 10mL liquid containing 1 of 9 different levels of human IgG calibrators
Manufacturer/Supplier:	ForteBio, Inc., 1360 Willow Road, Suite 205, Menlo Park, CA 94025-1516; Phone 650-322-1360, Fax 650-322-1370

### COMPOSITION / INFORMATION ON INGREDIENTS

Purified Human IgG in Phosphate Buffered Saline, pH 7.4, with 0.1% Bovine Serum Albumin, Tween 20 detergent (0.02%) and protease inhibitors. Sodium Azide (0.05%) used as preservative.

### HAZARDS IDENTIFICATION

Caution: The hazards of this material have not been thoroughly investigated. We recommend handling all materials with caution. The purified Human IgG was derived from source material which was tested and found negative for anti-HIV-1/HIV-2 (rDNA), negative for HIV-1 Antigen(s), negative for Anti-HCV, negative for Serological Test for Syphilis, and non-reactive for Hepatitis B Surface Antigen.

#### Potential Health Effects:

- Inhalation: Not determined
- Ingestion: Not determined
- Skin: Not determined
- Eyes: Not determined
- Chronic Exposure: Not determined
- Target Organs: Not determined

### FIRST AID MEASURES



1360 Willow Road, Suite 205  
Menlo Park, CA 94025-1516  
Phone 650-322-1360

## MATERIAL SAFETY DATA SHEET

For eye exposure rinse with water in an eyewash for at least 15 minutes with eyelids open. For skin exposure wash with soap and water; seek medical attention if irritation occurs. If swallowed, seek medical advice immediately.

### ACCIDENTAL RELEASE MEASURES

Collect and dispose of all waste in accordance with all applicable laws.

### HANDLING AND STORAGE

Store refrigerated (2 to 8°C).

### EXPOSURE CONTROL / PERSONAL PROTECTION

Wear gloves, protective clothing, and protective eyewear when handling the Calibrators and Negative Controls in this set. Follow safe laboratory practices.

ACGIH/OSHA Permissible Exposure Limit Data: Not determined

### PHYSICAL AND CHEMICAL PROPERTIES

Form: Liquid  
Odor: Not determined  
Solubility in Water: Not determined  
Specific Gravity: Not determined  
pH: 7.4  
Boiling Point: Not determined  
Melting Point: Not determined  
Flash Point: Not determined  
Vapor Pressure: Not determined

### STABILITY AND REACTIVITY

Thermal Decomposition: Not determined

Dangerous Reactions: Not determined



1360 Willow Road, Suite 205  
Menlo Park, CA 94025-1516  
Phone 650-322-1360

## MATERIAL SAFETY DATA SHEET

Dangerous Products of Decomposition: Not determined

### **TOXICOLOGICAL INFORMATION**

RTECS Number: None known  
Toxicity: Not determined  
Health Hazards: See above  
Potential Hazards: See above  
Carcinogenicity: Not determined.

### **ECOLOGICAL INFORMATION**

Dispose of according to all applicable laws.

### **DISPOSAL CONSIDERATIONS**

Consult local, state, and national regulations for disposal. Dispose of according to all applicable laws.

### **TRANSPORT INFORMATION**

Hazard Class: Not determined  
Identification Number: Not determined  
Packing Group: Not classified

Proper Shipping Name (Technical Name): Not classified

### **REGULATIONS**

US Toxic Substances Control Act (TSCA): Not listed  
US Other: Not applicable  
EEC EINECS Number: Not identified  
EEC Risk Statements: Not determined  
Other Country Regulations: None identified

### **OTHER INFORMATION**

This material is sold for Research Use Only. It is not intended for use in drug, household, agricultural nor cosmetics. Use of the product must be supervised by technically qualified individuals. The above information is correct to the best of our knowledge. Users should make their own independent decisions as regards completeness of information based on all sources possible. ForteBio, Inc., shall not be held liable for any damage resulting from handling or contact with the above product.