

24-Hour Emergency Number
(VelocityEHS) 1-800-255-3924 in U.S.

cleaning@alconox.com

alconox.com

- distributors
- technical information
- samples
- new developments
- white papers
- prices

TERGAZYME®

Enzyme-Active Powdered Detergent

- Concentrated to save you money
- Biodegradable and readily disposable
- Replaces corrosive acids and hazardous solvents
- Protease enzyme removes proteinaceous soils, tissue, blood and body fluids
- Free rinsing to provide reliable results and no interfering residues
- For FDA cleaning validations and food contact surface cleaning

Used to clean: Hospital instruments, dairy equipment, laboratory ware, reverse osmosis and ultrafiltration membranes and units, sampling apparatus, pharmaceutical and cosmetics manufacturing equipment, tubing, pipes, optical parts, process equipment, industrial parts, desalination plants, tanks and reactors.

Used to remove: Soils, grit, grime, blood, tissue, grease, fats, oils, proteinaceous soils, dairy proteins, particulates, solvents and bioreactor residue.

Surfaces cleaned: Corrosion inhibited formulation recommended for glass, metal, stainless steel, porcelain, ceramic, plastic, rubber and fiberglass. Can be used on soft metals such as copper, aluminum, zinc and magnesium if rinsed promptly. Corrosion testing may be advisable.

Cleaning method: Soak, brush, sponge, cloth, ultrasonic, flow through clean-in-place. Will foam—not for spray or machine use.

Directions: Make a fresh 1 –3% solution (3 – 8 tbs/gal, 1 1/4 – 4 oz/gal or 10 – 30 ml/L) in cold or warm water. If available, use warm water below 130°F (55°C). Clean by soak, circulate, wipe, or ultrasonic method. Follow manufacturers directions for filter membrane cleaning. Not for spray machines: will foam. RINSE THOROUGHLY—preferably with running water. For critical cleaning, do final or all rinsing in distilled, deionized, or purified water. For food contact surfaces, rinse with potable water. Used on a wide range of glass, ceramic, plastic, and metal

surfaces. Corrosion testing may be advisable. Material compatibility guidelines available on request.

Convenient Sizes	Alconox Cat. #
Case 9 x 4 lb (9 x 1.8 kg) jars	1304
25 lb (11 kg) box	1325
50 lb (22 kg) box	1350
100 lb (45 kg) drum	1301
300 lb (136 kg) drum	1303
1 lb makes 13 gal (1 kg makes 100L) cleaning solution	



TERGAZYME passes inhibitory residue test for water analysis and is FDA certified. TERGAZYME is available from leading laboratory, hospital, clinical and industrial suppliers. To find a distributor for Alconox Inc. detergents, use the Locate Dealers search tool at alconox.com. To request samples, use the Order Samples form at alconox.com. Or call Alconox Inc. at +914.948.4040, today.

TAZTB.5

24-Hour Emergency Number
(VelocityEHS) 1-800-255-3924 in U.S.

cleaning@alconox.com

alconox.com

- distributors
- technical information
- samples
- new developments
- white papers
- prices

Physical Data	Typical Value
pH of 1% Solution	9.5
Surface Tension 1% Sol'n (dyne/cm)	32
Conductivity 1% Sol'n (mS/cm at 21 degrees C)	6.3
Form	Powder
Color	White and cream colored flakes and brown specks
Solubility in Water	To 10% (w/w) at ambient temperature
Hard Water Effectiveness	Highly effective
Biodegradability	Yes, meets EC648/2004
Foam Tendency	High foaming
Shelf Life	42 months from end of month of manufacture
Fragrance Content	0%
Phosphate Content (as phosphorus)	17.4%
Flash Point (degrees F)	None
Organic Carbon (% calculated w/w)	11%

Chemical Description

Tergazyme consists primarily of a homogeneous blend of sodium linear alkylaryl sulfonate, phosphates, carbonates, and protease enzyme. Tergazyme contains anionic surfactants. The protease enzyme in Tergazyme is deactivated by 300 ppm hypochlorite at 85 degrees F in seconds; 3.5 ppm hypochlorite at 100 degrees F for 2 min; exposure to pH below 4 for 30 min at 140 degrees F; or by heating to 175 degrees F for 10 min.

Cleaning Validation and Residue Detection

To quantify trace levels of Tergazyme for cleaning validation purposes, use conductivity, UV, TOC, HPLC, or phosphorus detection. For details see the Cleaning Validation page in Support at alconox.com. For lab accreditation tests to show the detergent is rinsed away, see the Accreditation Residue Detection in the Support section of alconox.com.

Health and Safety Information

OSHA Hazardous Ingredients	None
RCRA Hazard Class	Nonhazardous
EPA Priority Pollutants	None
DOT Hazard Class	Nonhazardous
Flammability	Nonflammable
Latex Content	None in product or packaging.
Irritation	Cat. 2 skin, Cat. 2A eye irritant
Oral Toxicity	LD ₅₀ > 5000 mg/kg oral rat.
Inhalation Toxicity	Solution: No, Powder dust: irritant
VOC Content	0%
GMO Content	None
BSE/TSE Source	None
2007/68/EC Annex IIIa Allergens	None
REACH Art. 57 (SVHC)	None
RoHS	No RoHS hazardous ingredients.
Carcinogenicity	NTP = No IARC = No OSHA = No
Ingredients on TSCA Inventory	Yes
EC 907/2006	Contains enzymes; no optical brighteners, perfumes, allergenic fragrances, or preservative agents.

Precautions

Avoid contact with skin, eyes and clothing. Do not take internally. In case of contact immediately flush skin with plenty of water. For eyes, flush with plenty of water for at least 15 minutes and get medical attention immediately. A Safety Data Sheet is available at alconox.com. Contact Alconox Inc. for purchase specifications. Information presented is typical. Not to be taken as specifications. Typical data is not a specification.

While the information in this report should not be considered to be a product warranty, we urge you to investigate, test and verify the suitability of Alconox Inc. detergents for your specific application. We, of course, can not give permission to use, or recommend the use of, our detergents where they infringe patents. No representation or warranty is made as to the safety of products or materials mentioned under the Federal Food Additives Amendment of 1958.

TAZTB.5