

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

CITRANOX®

Acid Cleaner and Detergent

- Concentrated to save you money
- Phosphate-free, biodegradable and readily disposable
- Penetrating wetting power to save you time
- Free rinsing to give you reliable results and no interfering residues
- For FDA cleaning validations and food contact surface cleaning
- Contains acid ingredients for high performance removal of metal oxides, salts and inorganic residues combined with full detergent capability
- Powerful enough to be used for passivation of stainless steel

Used to clean: Dairy equipment, laboratory ware, sampling apparatus, lab stills, laser chillers, clean rooms, optical parts, electronic components, pharmaceutical and cosmetics manufacturing equipment, restaurant equipment, rethermalizers, steam tables, metal forgings, stampings and castings, industrial parts, pipes, tanks, boilers and reactors.

Used to remove: Scale, metal oxides, metal complexes, trace inorganics, milkstone, pasta and starch, soils, grit, buffing compound, grime, grease, fats, oils, particulates, deposits, amphoteric peptides, chemicals and solvents.

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, bronze, aluminum, and zinc if rinsed promptly. Corrosion testing may be advisable.

Cleaning method: Soak, brush, sponge, cloth, ultrasonic, immersion/recirculation cleaning in place. Will foam—not for spray or machine use.

Directions: Make a fresh 1–2% solution (3–5 tbsp per gal, 1 1/4–2 1/2 oz per gal, or 10–20 ml per L) in cold, warm or hot water. If available, use hot water. For difficult soils, raise water temperature and use more detergent. Clean by soak, circulate, wipe or ultrasonic method. Wear protective gloves and eyewear. 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.

For directions for passivation of stainless steel, search “Passivation” at alconox.com, or contact Alconox Inc. for the passivation technical bulletin.

Convenient Sizes	Alconox Cat. #
Case of 4 x 1 gal (4 x 3.8 L)	1801
5 gal (19 L) jerrycan	1805
15 gal (57 L) drum	1815
55 gal (208 L) drum	1855
1 gal makes 50 gal (1 L makes 50L) solution	



CITRANOX passes inhibitory residue test for water analysis and is FDA certified. CITRANOX 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.

CXTB.4

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 (as is)	2.5
Surface Tension 1% Sol'n (dyne/cm)	32
Conductivity 1% Sol'n (mS/cm at 21 degrees C)	0.7
Form	Liquid
Color	Pale yellow
Solubility in Water	Completely soluble in all proportions
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 and Dye Content	0%
Phosphate Content (as phosphorus)	0%
Flash Point (degrees F)	None
Organic Carbon (% calculated w/w)	17%
Density (lb/gal)	9.15

Chemical Description

Citranox contains a blend of organic acids, anionic and non-ionic surfactants and alkanolamines. Citranox is phosphate-free.

Cleaning Validation and Residue Detection

To quantify trace levels of Citranox for cleaning validation purposes, use conductivity, TOC, HPLC or anionic surfactant concentration. 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	Nonirritating
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 no enzymes, 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.

CXTB.4