

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

DETERGENT 8®

Low-Foaming Ion-Free Detergent

- Replace solvent cleaning
- Concentrated to save you money
- Biodegradable and readily disposable
- Penetrating wetting power to save you time
- Free rinsing to provide reliable results and no interfering residues
- Contains no phosphate, sulfur, chlorine, metal cation, halide, borate, silicate, carbonate, chlorocarbon, fluorocarbon or chelating ingredients

Used to clean: Circuit boards, electronic parts, phosphate sensitive labware, nuclear reactor cavities, delicate industrial parts and nuclear contaminated equipment.

Used to remove: Resins, rosins, oils, dirt, grime, particulates, residues, fluxes, chemicals and solvents.

Surfaces cleaned: Corrosion inhibited formulation recommended for glass, metal, stainless steel, ceramic, plastic, and circuit boards. Can be used, with prompt rinsing, on soft metals such as copper, aluminum, silver, and gold. Aluminum can darken and mixed metal galvanic reactions can occur. Corrosion testing is advisable.

Cleaning method: Soak, brush, sponge, cloth, ultrasonic, labware washer, spray washer, pressure washer and clean-in-place.

Directions: For manual, soak and ultrasonic use, make a 1 – 2% solution (3 – 5 Tbsp/gal, 1 1/4 – 2 1/2 oz/gal or 10 – 20 ml/L) in warm or hot water. An oily, milky solution is normal. Protective gloves and eyewear are recommended. For spray machines, add detergent at the hot water wash cycle to make a 2% solution (2 1/2 oz per gal, or 20 ml per liter). For difficult soils, raise water temperature and use more detergent.

RINSE THOROUGHLY—preferably with running water. For critical cleaning, do final or all rinsing in distilled, deionized or purified water. Material compatibility guidelines available on request.

Convenient Sizes Alconox Cat.

Case of 4 x 1 gal (4 x 3.8L)	1701
5 gal (19L) jerrycan	1705
15 gal (57L) drum	1715
55 gal (208L) drum	1755
1 gal makes 20–50 gal (1L makes 20–50L) solution	



DETERGENT 8 is FDA certified. DETERGENT 8 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.

D8TB.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 (1%)	11.0
Surface Tension 1% Sol'n (dyne/cm)	32
Conductivity 1% Sol'n (mS/cm at 21 degrees C)	0.7
Form	Liquid
Color	Clear
Solubility in Water	Completely soluble in all proportions
Hard Water Effectiveness	Highly effective
Biodegradability	Yes, meets EC648/2004
Foam Tendency	Low 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)	201
Organic Carbon (% calculated w/w)	35–40%
Density (lb/gal)	8.25

Chemical Description

Detergent 8 is an ion free, phosphate free, homogeneous blend of an alkanolamine, glycol ethers and an alkoxylated fatty alcohol. Contains no chelating agents, halides or conductive metal cations such as sodium. Detergent 8 contains non-ionic surfactants.

Cleaning Validation and Residue Detection

To quantify trace levels of Detergent 8 for cleaning validation purposes, Test a parameter of rinse water before and after rinsing the cleaned surface, or test the clean surface. No significant change in the parameter indicates no detectable residue. Parameters measured include spectrographic methods. For details, see the Cleaning Validation page in the Support section 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	UN1760, Corrosive Liquids, N.O.S. (Monoisopropanolamine Solution), 8, PG III
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	70% as is
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
Indredients 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.

D8TB.5