

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

Used to clean: Healthcare instruments, laboratory ware, sampling apparatus, plate glass, pharmaceutical and cosmetics manufacturing equipment, restaurant equipment, wine glasses, counters, floors, sidewalks and driveways, metal castings, forgings and stampings, industrial parts, tanks and reactors.

Used to remove: Soils, grit, grime, slime, grease, tars, resins, fats, oils, blood, tissue, inorganic residues, particulates, chemicals and solvents.

Surfaces cleaned: Corrosion inhibited formulation recommended for glass, metal, stainless steel, porcelain, ceramic, grout, 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: Labware washers, water-sterilizers, washer-sanitizers, cart washers, cage washers, spray washers, electrocleaning, ultrasonic cleaners, flow through clean-in-place. Manual cleaning if proper skin and eye protection is used.

Directions: Follow machine manufacturer's directions and use machine's powder detergent receptacles. In general, use 1 – 2% solution in hot water wash cycle (3 – 5 Tbsp/gal, 1 1/4 – 2 1/2 oz/gal or 10 – 20 ml/L). Most machines need 2 1/2 oz added to a typical 2 gal wash cycle. For manual, soak, circulate, or ultrasonic use, make a 1% solution in warm or hot water. For non-abrasive scrubbing,

make paste. For difficult soils, raise water temperature and use more detergent. Wear protective gloves and eyewear for manual cleaning. 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.

Convenient Sizes	Alconox Cat. #
Case 9 x 4 lb (9 x 1.8 kg) jars	1404
25 lb (11 kg) box	1425
50 lb (22 kg) box	1450
100 lb (45 kg) drum	1401
300 lb (136 kg) drum	1403
1 lb makes 26 gal (1 kg makes 200L) cleaning solution	



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

AJTB.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	11.0
Surface Tension 1% Sol'n (dyne/cm)	32
Conductivity 1% Sol'n (mS/cm at 21 degrees C)	8
Form	Powder
Color	White
Solubility in Water	Greater than 10% (w/w)
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 Content	0%
Phosphate Content (as phosphorus)	20%
Flash Point (degrees F)	None
Organic Carbon (% calculated w/w)	1.5%

Chemical Description

Alcojet is a blend of sodium silicate, a mixture of complex phosphate, concentrated organic chlorine, sodium carbonate, and an alkylalkoxylate. Alcojet is biodegradable and water soluble. Alcojet contains nonionic surfactants.

Cleaning Validation and Residue Detection

To quantify trace levels of Alcojet for cleaning validation purposes, use conductivity, 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 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.

AJT.B.5