

Pierce Sulfo-SANPAH, No-Weigh Format

MAN0017086

Rev. A.0

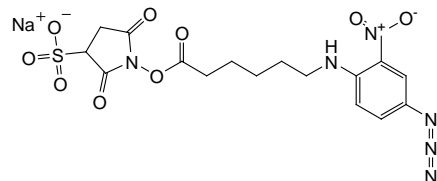
Pub. Part No. 2162637.0

A35395**Number****A35395****Description****Pierce Sulfo-SANPAH, No-Weigh Format** (sulfosuccinimidyl-6-[4'-azido-2'-nitrophenylamino]hexanoate), 10 × 1mg

Molecular Weight: 492.40

Spacer Arm: 18.2Å

Mass Added: 247.09



For Research Use Only. Not for use in diagnostic procedures.

Storage: Store Sulfo-SANPAH at -20°C. Store protected from moisture and light. Shipped on ice packs.**Introduction**

Thermo Scientific™ Pierce™ Sulfo-SANPAH is a heterobifunctional crosslinker that contains an amine-reactive *N*-hydroxysuccinimide (NHS) ester and a photoactivatable nitrophenyl azide. NHS esters react efficiently with primary amino groups (-NH₂) in pH 7-9 buffers to form stable amide bonds. The reaction results in the release of Sulfo-*N*-hydroxysuccinimide. When exposed to UV light, nitrophenyl azides form a nitrene group that can initiate addition reactions with double bonds, insertion into C-H and N-H sites, or subsequent ring expansion to react with a nucleophile (e.g., primary amines). Sulfo-SANPAH, supplied as a sodium salt, is useful for cell-surface protein crosslinking and possesses charged groups for water solubility to a concentration of 10mM.

Thermo Scientific™ No-Weigh™ products are specialty reagents provided in a pre-aliquoted format. The pre-weighed packaging prevents the loss of reagent reactivity and contamination over time by eliminating the repetitive opening and closing of the vial. The format enables use of a fresh vial of reagent each time, eliminating the hassle of weighing small amounts of reagents and reducing concerns over reagent stability.

Important Product Information

- Sulfo-SANPAH is moisture-sensitive. To avoid moisture condensation onto the product, equilibrate vial to room temperature before opening. Prepare immediately before use. The NHS-ester moiety readily hydrolyzes and becomes non-reactive; therefore, do not prepare stock solutions for storage. Discard any unused, reconstituted crosslinker.
- Hydrolysis of the NHS ester is a competing reaction and increases with increasing pH. Hydrolysis occurs more readily in dilute protein or peptide solutions. In concentrated protein solutions, the acylation reaction is favored.
- Use non-amine-containing buffers at pH 7-9 such as 20mM sodium phosphate, 0.15M NaCl (Product No. 28372); 20mM HEPES; 100mM carbonate/bicarbonate; or 50mM borate. Do not use buffers that contain Tris, glycine or sulfhydryls. Tris and glycine will compete with the intended reaction and thiols can reduce the azido group.
- For protein concentration greater than 5mg/mL, use a 10-fold molar excess of the crosslinker. For samples < 5mg/mL, use a 20- to 50-fold molar excess of the crosslinker. Use a final concentration of crosslinker at 0.1-10mM.

- Dissolve Sulfo-SANPAH in room-temperature water up to 10mM; solubility decreases with increasing salt concentration. Sulfo-SANPAH may also be dissolved in a dry water-miscible organic solvent such as DMSO or DMF. Add 200µL of buffer or solvent to a 1mg vial for a 10mM solution. Mix by repeated pipetting or replace cap and vortex. Higher molarity solutions can be obtained when DMSO or DMF is used rather than an aqueous buffer. The maximum useable volume of the vial is 800µL.
- For best results, react the NHS end of the crosslinkers (in the dark) first. After removing the hydrolyzed and non-reacted crosslinker by gel filtration or dialysis, the activated molecule can be coupled to a second molecule by photolysis.

Photolysis (Photoactivation) Information

- For maximum efficiency, use a shallow reaction vessel for photolysis. Irradiation efficiency decreases as the distance light must penetrate through the solution increases. Use a low protein-binding vessel for maximum sample recovery.
- For photolysis use a UV lamp that irradiates at 300-460nm (see **Note** below). High wattage lamps are more effective and require shorter exposure times than low wattage lamps. Suggestions for lamps include the Stratalinker 2400 (5 × 15 watt bulbs, emission at either 312nm or 365nm), mercury vapor lamps (180 watt, 350 watt, from 300 to 360nm), XeCl excimer laser (150mJ, 308nm) and UV Spectroline™ lamps (medium/long wavelength lamps). Using low wattage hand-held lamps, such as 6 watt lamps, will result in lower conjugation efficiencies.

Note: Avoid UV lamps that emit light at 254nm; this wavelength causes proteins to photodestruct. Filters that remove light at wavelengths below 300nm are ideal. Using a second filter that removes wavelengths above 370 nm could be beneficial but is not essential.

- Position the UV lamp 5-10cm from the reaction. For lamps >150 watts use a distance of 10cm. For lower powered lamps, use a distance of 5cm. Perform photolysis by placing the lamp above the reaction as the reaction vessel may impede irradiation by filtering some of the UV light.

Please visit our website for additional information relating to this product including the following items:

- Tech Tip #11: Light sources and conditions for photoactivation of aryl azide crosslinking reagents
- Tech Tip #43: Protein stability and storage
- Tech Tip #3: Determine reactivity of NHS-ester biotinylation and crosslinking reagents
- Tech Tip #5: Attach an antibody onto glass, silica or quartz surface

Related Thermo Scientific Products

21451	ANB-NOS (<i>N</i> -5-azido-2-nitrobenzoyloxysuccinimide), 50mg
22589	Sulfo-SANPAH (sulfosuccinimidyl-6-[4'-azido-2'-nitrophenylamino]hexanoate), 50mg
28372	BupH™ Phosphate Buffered Saline Pack, 40 packs
20290	DTT (Dithiothreitol), 5g
20291	No-Weigh™ DTT (Dithiothreitol), 48 × 7.7mg microtubes
35602	2-Mercaptoethanol
66382, 66807	Slide-A-Lyzer™ Dialysis Cassette Kits
66528	Slide-A-Lyzer™ Concentrating Solution, 200mL
66529	Slide-A-Lyzer Concentrating Solution, 500mL

General References

Gaudet, C., *et al.* (2003). Influence of type I collagen surface density on fibroblast spreading, motility, and contractility. *Biophys. J.* **85**:3329-35.
Uckun F.M., *et al.* (1995). Biotherapy of B-cell precursor leukemia by targeting genistein to CD19-associated tyrosine kinases. *Science* **267**:886-91.

Products are warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation"). No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

NO OTHER WARRANTIES, EXPRESS OR IMPLIED, ARE GRANTED, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR NON INFRINGEMENT. BUYER'S EXCLUSIVE REMEDY FOR NON-CONFORMING PRODUCTS DURING THE WARRANTY PERIOD IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR PRODUCTS AS THE RESULT OF (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, (II) MISUSE, FAULT OR NEGLIGENCE OF OR BY BUYER, (III) USE OF THE PRODUCTS IN A MANNER FOR WHICH THEY WERE NOT DESIGNED, OR (IV) IMPROPER STORAGE AND HANDLING OF THE PRODUCTS.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at thermofisher.com. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2017 Thermo Fisher Scientific Inc. All rights reserved. Spectroline is a trademark of Spectronics Corp. All (other) trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.