

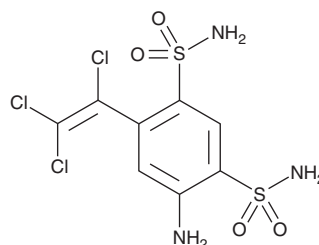
# PRODUCT INFORMATION



## Clorsulon

Item No. 30288

**CAS Registry No.:** 60200-06-8  
**Formal Name:** 4-amino-6-(1,2,2-trichloroethenyl)-1,3-benzenedisulfonamide  
**Synonyms:** L-631529, MK 401  
**MF:** C<sub>8</sub>H<sub>8</sub>Cl<sub>3</sub>N<sub>3</sub>O<sub>4</sub>S<sub>2</sub>  
**FW:** 380.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 227, 268, 325 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Clorsulon is supplied as a solid. A stock solution may be made by dissolving the clorsulon in the solvent of choice, which should be purged with an inert gas. Clorsulon is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of clorsulon in these solvents is approximately 10 mg/ml in ethanol and DMSO and approximately 15 mg/ml in DMF.

Clorsulon is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, clorsulon should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Clorsulon has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Clorsulon is a sulfonamide antiparasitic agent.<sup>1</sup> It induces swelling and blebbing and reduces motility of *F. hepatica* mature flukes when used at a concentration of 10 µg/ml. *In vivo*, clorsulon (3.5-15 mg/kg) reduces the number of flukes in a goat model of *F. hepatica* infection.<sup>2</sup> It completely eradicates naturally acquired *F. hepatica* adult flukes in sheep and cattle when administered at a dose of 7 mg/kg.<sup>3</sup> Formulations containing clorsulon have been used in the treatment of liver flukes in sheep and cattle.

### References

1. Meaney, M., Fairweather, I., Brennan, G.P., *et al.* *Fasciola hepatica*: Effects of the fasciolicide clorsulon in vitro and in vivo on the tegumental surface, and a comparison of the effects on young- and old-mature flukes. *Parasitol. Res.* **91**(3), 238-250 (2003).
2. Sundlof, S.F., Bliss, E.L., Greiner, E.C., *et al.* Efficacy of clorsulon for the treatment of experimentally induced infections of *Fasciola hepatica* in goats. *Am. J. Vet. Res.* **52**(1), 111-114 (1991).
3. Zimmerman, G.L., Wallace, D.H., Schons, D.J., *et al.* Efficacy of clorsulon against mature naturally acquired *Fasciola hepatica* infections in cattle and sheep. *Am. J. Vet. Res.* **47**(8), 1665-1667 (1986).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/08/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)