

PRODUCT DATA SHEET



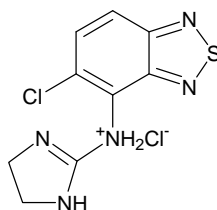
PRODUCT: Tizanidine·HCl
5-Chloro-4-(2-imidazolin-2-yl-amino)-
2,1,3-benzothiadiazol·HCl

CAS NO: 51322-75-9

CATALOG NO.: AR-103

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid

MOLECULAR FORMULA: C₉H₈ClN₅S·HCl

MOLECULAR WEIGHT: 290.2

PURITY: 99%

SOLUBILITY: Soluble in water (20 mg/ml)

STORAGE: Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.

APPLICATION NOTES: α_2 -adrenergic receptor agonist.^{1,2} Displays affinity for the I₁-imidazoline receptor.³ Spasmolytic activity.⁴

- REFERENCES:
1. T.Kameyama *et al. Eur.J.Pharmacol.* 1986 **125** 257
 2. T.Nabeshima *et al. Neuropharmacol.* 1987 **26** 1453
 3. J.E.Piletz *et al. J.Pharmacol.Exp.Ther.* 1996 **26** 462
 4. D.M.Coward *Neurology* 1994 **44** S6

The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.