

## NSC-207895/XI-006

### 2,1,3-Benzoxadiazole,4-(4-methyl-1-piperazinyl)-7-nitro-,3-oxide

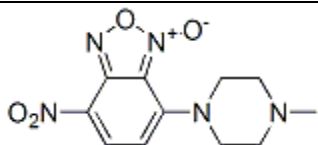
Cat. # SI9825

**Background:** A small molecule inhibitor that has been found to serve as an inhibitor of MDMX. NSC-207895 has shown capability of inducing apoptosis in MCF-7 cells.<sup>1</sup>

**Application:** Inhibits MDMX expression resulting in increased p53 activity and DNA damage

#### Product Information:

<b>CAS No.</b>	58131-57-0
<b>Purity:</b>	>99% by HPLC
<b>Molecular Weight:</b>	279.25 Da
<b>Physical State:</b>	Powder
<b>Quantity:</b>	5mg
<b>Solubility:</b>	DMSO (0.4 mg/ml); Ethanol (<1mg/ml); Water (<1mg/ml)
<b>Storage:</b>	Store desiccated as supplied at -20°C for 2 years.



#### Formula:

C<sub>11</sub>H<sub>13</sub>N<sub>5</sub>O<sub>4</sub>

#### References

1. Wang H, et al. (2011). "A small-molecule inhibitor of MDMX activates p53 and induces apoptosis." *Mol Cancer Ther* 10(1): 69-79.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2014 LifeSensors, Inc. All Rights Reserved