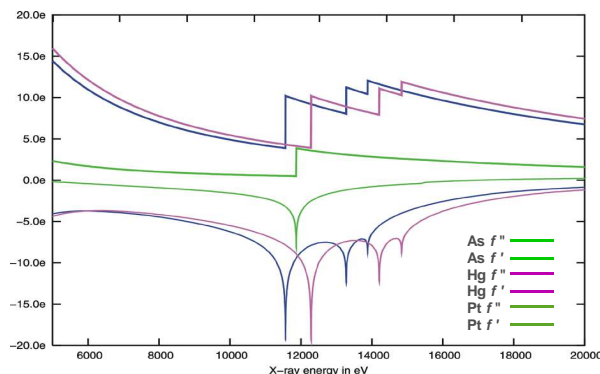
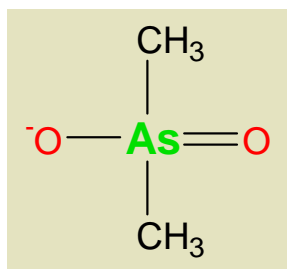
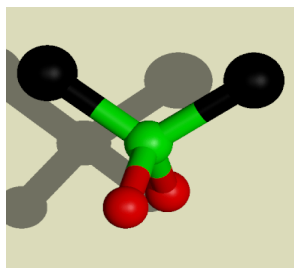


# Cacodylate is very toxic



- **can interfere with your MAD experiment**

The K absorption edge of As is at an energy of 11.8667 keV, close to the LIII absorption edges of Hg (12.2839 keV) and Pt (11.5637 keV).



- **can distort the active site of enzymes [1]**

since it is able to react with free sulfhydryl groups and thus can covalently bind to cysteine sulfur. Your protein conformation may be seriously affected [2]!

## HEALTH HAZARD INFORMATION

**May be fatal if swallowed. Known carcinogen in humans.**

Harmful if inhaled, may be harmful by skin contact. Long-term exposure may lead to kidney and liver damage. Eye and skin irritant.

**Cacodylate (pK<sub>a</sub> 6.3) is substituted by MES (pK<sub>a</sub> 6.2) in all JBScreens!**

[1] Goldgur *et al.* (1998) Three new structures of the core domain of HIV-1 integrase: An active site that binds magnesium. *Proc. Natl. Acad. Sci.* **95**:9150.

[2] Noel *et al.* (1993) The 2.2 Å crystal structure of transducin-α complexed with GTP γS. *Nature* **366**:654.