

## Cholinesterase References

### Reference

1. Huang, X., Lee, B. W., Johnson, G. L., Naleway, J., Guzikowski, A., Dai, W., Darzynkiewicz, Z. 2005. Novel assay utilizing fluorochrome-tagged physostigmine (Ph-F) to in situ detect active acetylcholinesterase (AChE) induced during apoptosis. *Cell Cycle* 4:1, 140-147.