



## Isotype composition of isolated tubulins

### Beta-tubulin Isotype composition of Mammalian Brain and Cancer Cell Line Sources.

Source	Tubulin isotype				Reference
	$\beta$ I/ $\alpha$	$\beta$ II/ $\alpha$	$\beta$ III/ $\alpha$	$\beta$ IV/ $\alpha$	
Bovine brain	3%	58%	25%	13%	Banjeree and Luduena 1992
HeLa S3 suspension ATCC CCL-2.2	80%	0%	0%	20%	Newton et al. 2002
HeLa S3 suspension ATCC CCL-2.2	90%	0%	0%	10%	Davis et al. 2009
HeLa attached ATCC CCL-2	90%	0%	0%	10%	Davis et al. 2009
MCF-7 cell line ATCC HTB-22	39%	0%	3%	58%	Hiser et al. 2006
MCF-7 cell line	55%	0%	6%	39%	Davis et al. 2009



## Isotype composition of isolated tubulins

### References

Banerjee, A., and Luduena, R. F. 1992. Kinetics of colchicine binding to purified  $\beta$ -tubulin isotypes from bovine brain. *J. Biol. Chem.* 267, 13335–13339.

Davis A, Martinez S, Nelson D. and Middleton K. 2009. A Tubulin Polymerization Microassay Used to Compare Ligand Efficacy. *Meth. Cell Biol.* 95, Ch 18, p.327-347.

Hiser, L., Aggarwal, A., Young, R., Frankfurter, A., Spano, A., Correia, J. J., and Lobert, S. 2006. Comparison of beta-tubulin mRNA and protein levels in 12 human cancer cell lines. *Cell Motil. Cytoskeleton*, 63, 41–52.

Newton, C. N., DeLuca, J. G., Himes, R. H., Miller, H. P., Jordan, M. A., and Wilson, L. 2002. Intrinsically slow dynamic instability of HeLa cell microtubules in vitro. *J. Biol. Chem.* 277, 42456–42462.