

Montana Molecular's cAMP sensor in HEK-293 cells

