

Multiplex Ca<sup>2+</sup> and cAMP in HEK-293 Cells

