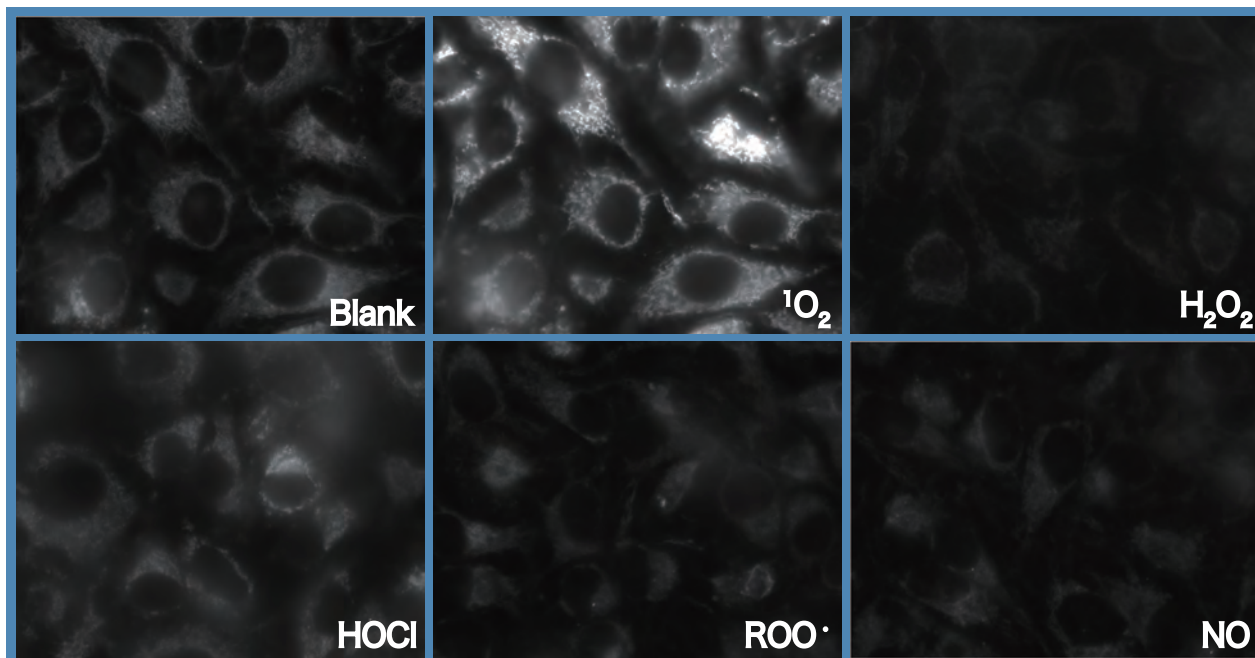
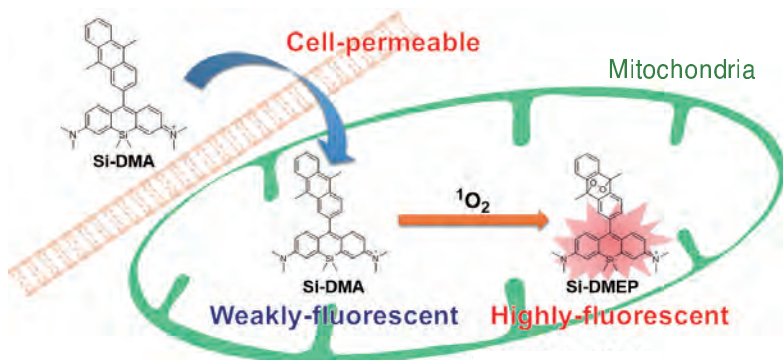


Si-DMA_{for} Mitochondrial Singlet Oxygen Imaging

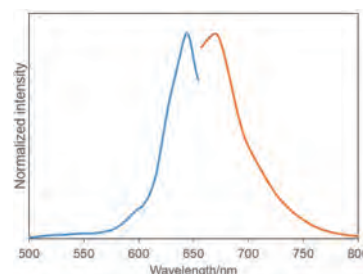
DOJINDO MOLECULAR TECHNOLOGIES, INC.



Selectivity of Si-DMA toward various ROS in living cells



Singlet oxygen detection mechanism by Si-DMA



Excitation and emission spectra of Si-DMEP



Si-DMA is available at **Dojindo**

Poster Session

2016. 4.7 Thr

16:20-17:50

No. 4P09

on the 3rd floor of the venue

URL: <http://www.dojindo.co.jp/>

E-mail: info@dojindo.co.jp