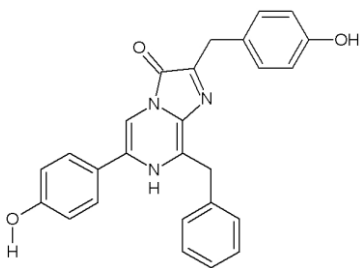
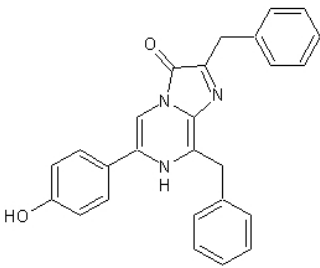
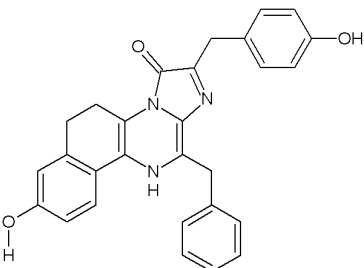
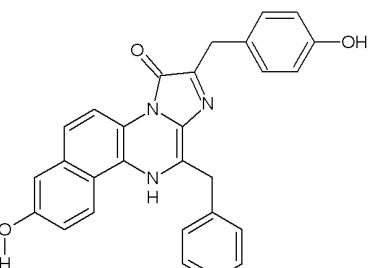
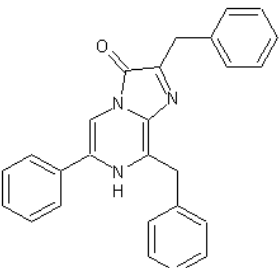
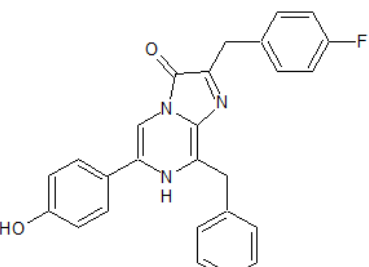
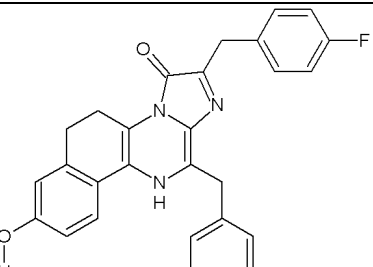
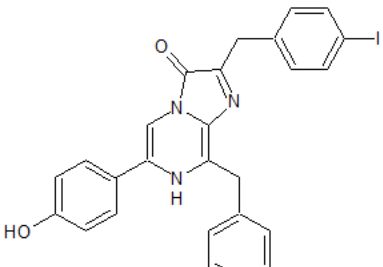


Compound	Cat. #	Structure	Abbrev.	M.W.	total light intensity & emission max with:							
					RmLuc		GLuc		NLuc		Aequorin	
					% light	Em _{max}	% light	Em _{max}	% light	Em _{max}	% light	Em _{max}
Coelenterazine (native)	303		CTZ	423.5	100	475	100	480	100	457	100	465
h-Coelenterazine	301		h-CTZ	407.5	41	475	0	n/a	1060	455	220	464
e-Coelenterazine	355		e-CTZ	449.5	137	418/ 475	0	n/a	UNK	UNK	50	405/ 465
v-Coelenterazine	370		v-CTZ	447.5	73	520	0	n/a	UNK	UNK	3.5	UNK
Coelenterazine 400a	340		CTZ-400a	391.5		400	0	n/a	770	453	UNK	UNK
Coelenterazine-F	345		CTZ-F	425.5	58/ 122	472	0	n/a	920	453	270	473
e Coelenterazine-F	346		eCTZ-F	451.5	130	413/ 475	6	n/a	UNK	UNK	UNK	UNK
Coelenterazine-I	347		CTZ-I	533.4	5	UNK	18	UNK	UNK	UNK	UNK	UNK

*information based on following data: 1) Inouye S, Shimomura O. Biochem Biophys Res Commun. 1997 Apr 17;233(2):349-53.
2) Inouye S, et al. Biochem Biophys Res Commun. 2013 Jul 19;437(1):23-8.
3) unpublished data by Prolume