

Transcreener Assay Setup Guide on the BMG LABTECH VANTASTAR™ Microplate Reader

Transcreener FI

| | |
|--------------------------|---|
| Filters | |
| Excitation | 580-10 nm |
| Dichroic | 601 nm |
| Emission | 620-10 nm |
| Gain | obtained through adjustment on tracer only control, here ~ 1770 |
| Focal height | 12,2, focus adjustment performed on tracer only control |
| Settling time | 0.1 s |
| Number of flashes | 100 |

Transcreener FP

| | |
|--------------------------|--|
| Filters | |
| Excitation | 590-50 nm |
| Dichroic | LP 639 |
| Emission 1 | 675-50 |
| Gain | Was set to achieve 20 mP for a well containing the free tracer, here ~1300 |
| Focal height | 11.6, focus adjustment performed on positive control |
| Settling time | 0.1 s |
| Number of flashes | 100 |

Transcreener TR-FRET

| | |
|--------------------------|---|
| Filters | |
| Excitation | 337 nm, Ex TR |
| Dichroic | Included, LP TR |
| Emission 1 | 620-10 nm |
| Emission 2 | 665-10 nm |
| Gain | obtained through adjustment on S1 (100% ATP conversion) for 620nm channel (donor), gain value transferred to 665nm channel (acceptor), here ~1800 |
| Focal height | 11.2, focus adjustment performed on positive control |
| Settling time | 0.3 s |
| Number of flashes | 100 |
| Integration Start | 60 μ s |
| Integration time | 400 μ s |