

Datasheet of GLuc-ONTM PKC Transcriptional Response Element (TRE)  
Clone (TR106)

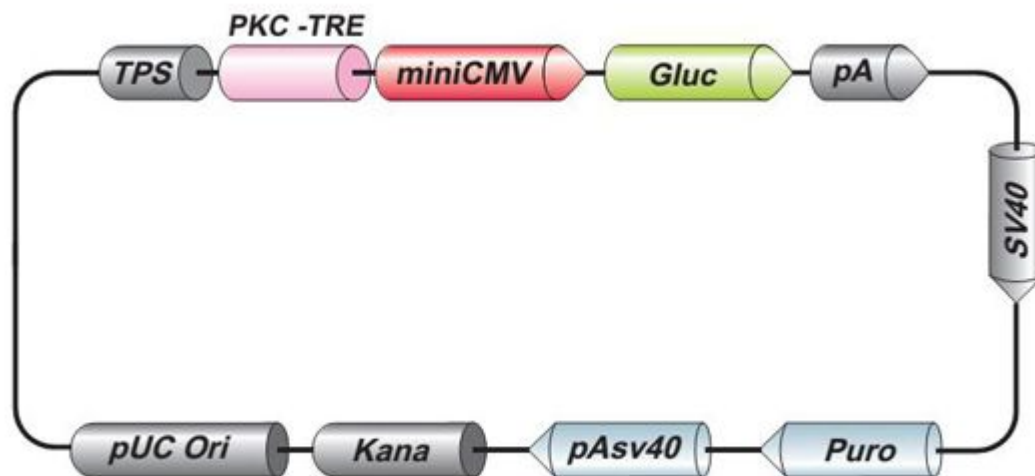


Figure 1. GLuc-ONTM PKC Transcriptional Response Element (TRE) Clone(TR106)

**Catalog:**

TR106

**Description:**

The PKC transcriptional response element (TRE) clone assays the status of the PKC/MAPK signaling pathway through the activity of the transcription factor PKC. The PKC transcriptional TRE clone contains a minimal promoter and tandem repeats of the PKC transcriptional response element upstream of a secreted Gaussia luciferase reporter gene. The number of response elements, as well as the intervening sequence between the elements has been optimized to maximize the sensitivity and signal to noise ratio.

**Delivery:**

500ng/μL × 40 μL

**Storage Conditions:**

-20°C

**QC:**

- . Functional validation in common cell line
- . Full sequence verification
- . Restriction Digestion

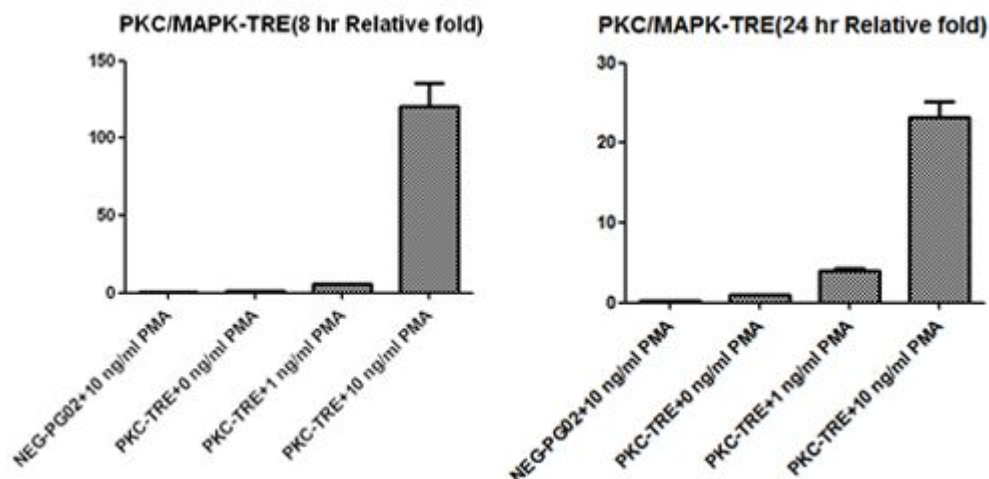


Figure 2. Validation data of GLuc-ON™ PKC Transcriptional Response Element (TRE) Clone(TR106).

## Restriction Enzyme Information for TR106

Table 1. Restriction Enzymes That Do Not Cleave TR106

|       |         |        |       |       |       |       |       |
|-------|---------|--------|-------|-------|-------|-------|-------|
| AfeI  | AgeI    | AscI   | BaeI  | BaeI  | BbvCI | BlpI  | BsaAI |
| BstXI | BstZ17I | Bsu36I | BtrI  | EcoRV | FspI  | MluI  | NdeI  |
| PacI  | PmeI    | PmlI   | PpuMI | PshAI | PvuII | SanDI | SapI  |
| SbfI  | SgrAI   | SnaBI  | SpeI  | SrfI  | Swal  |       |       |

Table 2. Restriction Enzymes That Cleave TR106 Once

|         |      |         |      |         |      |         |      |        |      |          |      |          |      |          |      |
|---------|------|---------|------|---------|------|---------|------|--------|------|----------|------|----------|------|----------|------|
| AarI    | 619  | AasI    | 4526 | AccI    | 1190 | AccIII  | 2166 | AcII   | 209  | AcuI     | 4086 | Adel     | 2624 | AflII    | 343  |
| AflIII  | 4628 | AhdI    | 1155 | Alol    | 880  | Aor13HI | 2166 | AseI   | 3709 | AsiSI    | 3510 | AspA2I   | 1976 | AspEI    | 1155 |
| AspI    | 2095 | AssI    | 554  | AsuII   | 458  | AvrII   | 1976 | BamHI  | 840  | BauI     | 4455 | BciVI    | 4430 | BclI     | 2027 |
| BfrI    | 343  | BfuI    | 4430 | BlnI    | 1976 | BmcAI   | 554  | BmeRI  | 1155 | Bpu14I   | 458  | Bsa29I   | 547  | BsaBI    | 883  |
| BsaI    | 2527 | Bse8I   | 883  | BseAI   | 2166 | BseCI   | 547  | BseJI  | 883  | BsePI    | 2508 | BseRI    | 1972 | BshVI    | 547  |
| BsiWI   | 2109 | Bso31I  | 2527 | Bsp119I | 458  | Bsp13I  | 2166 | BspDI  | 547  | BspEI    | 2166 | BspHI    | 3034 | BspT104I | 458  |
| BspTI   | 343  | BssHII  | 2508 | BssSI   | 4455 | Bst2BI  | 4455 | BstAFI | 343  | BstBI    | 458  | BstEII   | 2187 | BstENI   | 3422 |
| BstPI   | 2187 | Bsu15I  | 547  | Bsul    | 4430 | CciI    | 3034 | CciNI  | 1412 | Cfr9I    | 1997 | ClaI     | 547  | CpoI     | 2169 |
| CspI    | 2169 | DraI    | 2934 | DraII   | 2624 | DrdI    | 4526 | DriI   | 1155 | DseDI    | 4526 | Eam1105I | 1155 | Ecl136II | 672  |
| Eco31I  | 2527 | Eco53kI | 672  | Eco57I  | 4086 | Eco91I  | 2187 | EcoCRI | 672  | EcoNI    | 3422 | EcoO65I  | 2187 | EcoRI    | 483  |
| FbaI    | 2027 | FblI    | 1190 | FseI    | 2732 | HindIII | 764  | HpaI   | 2873 | Kpn2I    | 2166 | KroI     | 2728 | Ksp22I   | 2027 |
| KspAI   | 2873 | MfeI    | 2882 | MroI    | 2166 | MroNI   | 2728 | MspCI  | 343  | MunI     | 2882 | NaeI     | 2730 | NgoMIV   | 2728 |
| NotI    | 1412 | NspV    | 458  | PaeR7I  | 1418 | PagI    | 3034 | PasI   | 2102 | Paul     | 2508 | PciI     | 4628 | PdiI     | 2730 |
| Pfi23II | 2109 | PfiFI   | 2095 | PfoI    | 2528 | PscI    | 4628 | PshBI  | 3709 | Psp124BI | 674  | Psp1406I | 209  | PspEI    | 2187 |
| PspLI   | 2109 | PspXI   | 1418 | PstI    | 2975 | PsyI    | 2095 | PteI   | 2508 | RcaI     | 3034 | RgaI     | 3510 | RigI     | 2732 |
| Rsr2I   | 2169 | RsrII   | 2169 | SacI    | 674  | Sall    | 1189 | Scal   | 554  | SfaAI    | 3510 | Sfr274I  | 1418 | SfuI     | 458  |
| SgfI    | 3510 | Slal    | 1418 | SmaI    | 1999 | SspI    | 3435 | SstI   | 674  | StrI     | 1418 | TspMI    | 1997 | Tth111I  | 2095 |
| Vha464I | 343  | VspI    | 3709 | XagI    | 3422 | XbaI    | 1430 | XcmI   | 1651 | XhoI     | 1418 | XmaI     | 1997 | XmaJI    | 1976 |
| XmiI    | 1190 | Zrml    | 554  |         |      |         |      |        |      |          |      |          |      |          |      |

Table 3. Restriction Enzymes That Cleave TR106 Twice

|         |      |         |      |         |      |         |      |          |      |          |      |          |      |          |      |
|---------|------|---------|------|---------|------|---------|------|----------|------|----------|------|----------|------|----------|------|
| Acc36I  | 619  | Acc36I  | 1325 | Acc65I  | 663  | Acc65I  | 2709 | AccB7I   | 479  | AccB7I   | 3773 | Ajul     | 462  | Ajul     | 3538 |
| Alw44I  | 1095 | Alw44I  | 4314 | AlwNI   | 1067 | AlwNI   | 4219 | Apal     | 356  | Apal     | 3936 | ApaLI    | 1095 | ApaLI    | 4314 |
| Asp700I | 1114 | Asp700I | 1438 | Asp718I | 663  | Asp718I | 2709 | AsuNHI   | 44   | AsuNHI   | 676  | Ball     | 946  | Ball     | 2428 |
| BasI    | 479  | BasI    | 3773 | BbsI    | 134  | BbsI    | 929  | Bcgl     | 71   | Bcgl     | 2494 | BfuAI    | 619  | BfuAI    | 1325 |
| BglII   | 475  | BglII   | 745  | BmtI    | 48   | BmtI    | 680  | Bpil     | 134  | Bpil     | 929  | BpII     | 1911 | BpII     | 1943 |
| Bpu10I  | 668  | Bpu10I  | 3527 | BpuAI   | 134  | BpuAI   | 929  | BpvUI    | 1187 | BpvUI    | 3510 | BsaXI    | 878  | BsaXI    | 2431 |
| BsmBI   | 2445 | BsmBI   | 3526 | Bsp120I | 352  | Bsp120I | 3932 | Bsp1407I | 440  | Bsp1407I | 1552 | Bsp19I   | 851  | Bsp19I   | 1883 |
| Bsp68I  | 958  | Bsp68I  | 3167 | BspMI   | 619  | BspMI   | 1325 | BspOI    | 48   | BspOI    | 680  | BsrGI    | 440  | BsrGI    | 1552 |
| BstAUI  | 440  | BstAUI  | 1552 | BstV2I  | 134  | BstV2I  | 929  | BtgZI    | 870  | BtgZI    | 1227 | BtuMI    | 958  | BtuMI    | 3167 |
| BveI    | 619  | BveI    | 1325 | Cail    | 1067 | Cail    | 4219 | Cfr42I   | 984  | Cfr42I   | 2267 | Csil     | 1743 | Csil     | 2620 |
| DinI    | 2262 | DinI    | 2393 | Drall   | 353  | Drall   | 3933 | Eco147I  | 1975 | Eco147I  | 2384 | EcoO109I | 353  | EcoO109I | 3933 |
| EgeI    | 2262 | EgeI    | 2393 | Ehel    | 2262 | Ehel    | 2393 | Esp3I    | 2445 | Esp3I    | 3526 | HincII   | 1191 | HincII   | 2873 |
| HindII  | 1191 | HindII  | 2873 | KasI    | 2260 | KasI    | 2391 | KpnI     | 667  | KpnI     | 2713 | KspI     | 984  | KspI     | 2267 |
| MabI    | 1743 | MabI    | 2620 | MIsI    | 946  | MIsI    | 2428 | MluNI    | 946  | MluNI    | 2428 | Mly113I  | 2261 | Mly113I  | 2392 |
| MroXI   | 1114 | MroXI   | 1438 | MscI    | 946  | MscI    | 2428 | MslI     | 524  | MslI     | 889  | Msp20I   | 946  | Msp20I   | 2428 |
| MvrI    | 1187 | MvrI    | 3510 | NarI    | 2261 | NarI    | 2392 | NcoI     | 851  | NcoI     | 1883 | NheI     | 44   | NheI     | 676  |
| NruI    | 958  | NruI    | 3167 | PceI    | 1975 | PceI    | 2384 | Pdml     | 1114 | Pdml     | 1438 | PfiMI    | 479  | PfiMI    | 3773 |
| Ple19I  | 1187 | Ple19I  | 3510 | PluTI   | 2264 | PluTI   | 2395 | PspOMI   | 352  | PspOMI   | 3932 | PstNI    | 1067 | PstNI    | 4219 |
| PvuI    | 1187 | PvuI    | 3510 | RruI    | 958  | RruI    | 3167 | RseI     | 524  | RseI     | 889  | SacII    | 984  | SacII    | 2267 |
| SexAI   | 1743 | SexAI   | 2620 | SfiI    | 657  | SfiI    | 758  | SfoI     | 2262 | SfoI     | 2393 | Sfr303I  | 984  | Sfr303I  | 2267 |
| SgrBI   | 984  | SgrBI   | 2267 | SmiMI   | 524  | SmiMI   | 889  | SseBI    | 1975 | SseBI    | 2384 | SspDI    | 2260 | SspDI    | 2391 |
| StuI    | 1975 | StuI    | 2384 | Van91I  | 479  | Van91I  | 3773 | VneI     | 1095 | VneI     | 4314 | XmnI     | 1114 | XmnI     | 1438 |

Table 4. Restriction Enzymes That Cleave TR106 Three Times

|          |      |          |      |          |      |          |      |          |      |        |      |        |      |          |      |
|----------|------|----------|------|----------|------|----------|------|----------|------|--------|------|--------|------|----------|------|
| AanI     | 427  | AanI     | 1569 | AanI     | 2853 | AatII    | 66   | AatII    | 2099 | AatII  | 2597 | Avall  | 2169 | Avall    | 2279 |
| Avall    | 2611 | BfmI     | 2971 | BfmI     | 4172 | BfmI     | 4363 | Bme18I   | 2169 | Bme18I | 2279 | Bme18I | 2611 | Bmrl     | 141  |
| Bmrl     | 611  | Bmrl     | 1855 | Bmul     | 141  | Bmul     | 611  | Bmul     | 1855 | Bpml   | 1033 | Bpml   | 1252 | Bpml     | 2551 |
| Bse3DI   | 421  | Bse3DI   | 1577 | Bse3DI   | 3008 | BseMI    | 421  | BseMI    | 1577 | BseMI  | 3008 | BsrDI  | 421  | BsrDI    | 1577 |
| BsrDI    | 3008 | Bst6I    | 329  | Bst6I    | 2212 | Bst6I    | 3323 | BstAPI   | 1721 | BstAPI | 1793 | BstAPI | 2624 | BstSFI   | 2971 |
| BstSFI   | 4172 | BstSFI   | 4363 | Eam1104I | 329  | Eam1104I | 2212 | Eam1104I | 3323 | EarI   | 329  | EarI   | 2212 | EarI     | 3323 |
| Eco47I   | 2169 | Eco47I   | 2279 | Eco47I   | 2611 | Gsul     | 1033 | Gsul     | 1252 | Gsul   | 2551 | Psil   | 427  | Psil     | 1569 |
| Psil     | 2853 | Sfcl     | 2971 | Sfcl     | 4172 | Sfcl     | 4363 | TaqII    | 2071 | TaqII  | 2695 | TaqII  | 3764 | VpaK11BI | 2169 |
| VpaK11BI | 2279 | VpaK11BI | 2611 | Zral     | 64   | Zral     | 2097 | Zral     | 2595 |        |      |        |      |          |      |