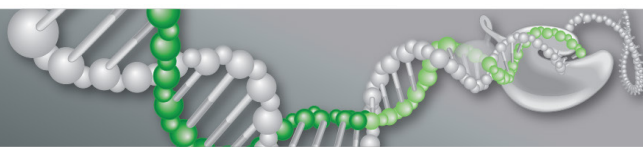


PrimePCR™ Pricing

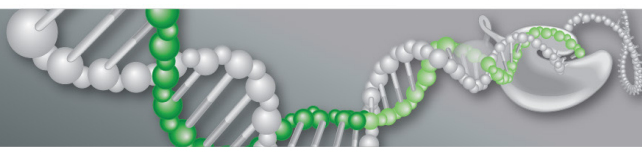


アッセイおよびコントロール

プライマー アッセイ (SYBR Green assay) 価格

| Product | Catalog # | Description | List Price (JPY) |
|---|---------------------|--|------------------|
| mRNA Primer Assays (desalted) (Wet-lab validated — human, mouse, and rat) | 10025636 | Primer assay desalted, 200 reactions | ¥ 16,000 |
| | 10025637 | Primer assay desalted, 1,000 reactions | ¥ 48,000 |
| | 10025638 | Primer assay desalted, 2,500 reactions | ¥ 80,000 |
| mRNA Primer Assays (HPLC) (Wet-lab validated — human, mouse, and rat) | 10025639 | Primer assay HPLC, 200 reactions | ¥ 26,000 |
| | 10025640 | Primer assay HPLC, 1,000 reactions | ¥ 78,000 |
| | 10025641 | Primer assay HPLC, 2,500 reactions | ¥ 130,000 |
| mRNA Primer Assays (desalted) (Expertly designed — Arabidopsis, chicken, cow, dog, pig, rabbit, rhesus monkey, yeast, zebrafish) | 10042976 | Primer assay desalted, 200 reactions | ¥ 12,000 |
| | 10042977 | Primer assay desalted, 1,000 reactions | ¥ 36,000 |
| | 10042978 | Primer assay desalted, 2,500 reactions | ¥ 60,000 |
| mRNA Primer Assays (HPLC) (Expertly designed — Arabidopsis, chicken, cow, dog, pig, rabbit, rhesus monkey, yeast, zebrafish) | 10042979 | Primer assay HPLC, 200 reactions | ¥ 24,000 |
| | 10042980 | Primer assay HPLC, 1,000 reactions | ¥ 72,000 |
| | 10042981 | Primer assay HPLC, 2,500 reactions | ¥ 120,000 |
| Custom Primer Pairs (desalted) | 10025220 | Custom primer pair desalted, 200 reactions | ¥ 6,000 |
| | 10025221 | Custom primer pair desalted, 1,000 reactions | ¥ 16,000 |
| | 10025222 | Custom primer pair desalted, 2,500 reactions | ¥ 26,000 |
| Custom Primer Pairs (HPLC) | 10025223 | Custom primer pair HPLC, 200 reactions | ¥ 22,000 |
| | 10025224 | Custom primer pair HPLC, 1,000 reactions | ¥ 66,000 |
| | 10025225 | Custom primer pair HPLC, 2,500 reactions | ¥ 110,000 |
| PreAmp Assays | 10041594 | PreAmp custom assay, 400 reactions | ¥ 4,000 |
| | 10041595 | PreAmp SYBR® Green assay, 400 reactions | ¥ 16,000 |
| | 10041596 | PreAmp probe assay, 400 reactions | ¥ 16,000 |
| lncRNA Primer Assays (desalted) (Expertly designed — human) | 12004175 | lncRNA Primer assay desalted, 200 reactions | ¥ 16,000 |
| | 12004176 | lncRNA Primer assay desalted, 1,000 reactions | ¥ 48,000 |
| | 12004177 | lncRNA Primer assay desalted, 2,500 reactions | ¥ 80,000 |
| lncRNA Primer Assays (HPLC) (Expertly designed — human) | 12004178 | lncRNA Primer assay HPLC, 200 reactions | ¥ 26,000 |
| | 12004179 | lncRNA Primer assay HPLC, 1,000 reactions | ¥ 78,000 |
| | 12004180 | lncRNA Primer assay HPLC, 2,500 reactions | ¥ 130,000 |
| lncRNA PreAmp Assays | 12004197 | lncRNA PreAmp SYBR® Green assay, 400 reactions | ¥15,000 |
| | 12004198 | lncRNA PreAmp probe assay, 400 reactions | ¥15,000 |
| PreAmp Pool for lncRNA Arrays | 対応するアレイの種類によって異なります | lncRNA PreAmp pool, 25 reactions | ¥ 32,000 |

PrimePCR™ Pricing



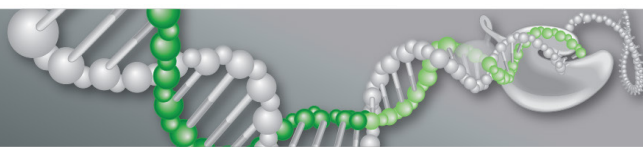
アッセイおよびコントロール

プローブ アッセイ価格

| Product | Catalog # | Description | List Price (JPY) |
|--|---|--|------------------|
| mRNA qPCR Probe Assays (Wet-lab validated — human, mouse, and rat) | 12001950 | qPCR FAM probe assay, 200 reactions | ¥ 24,000 |
| | 10031225 | qPCR FAM probe assay, 500 reactions | ¥ 44,000 |
| | 10031226 | qPCR FAM probe assay, 1,000 reactions | ¥ 62,000 |
| | 10031227 | qPCR FAM probe assay, 2,500 reactions | ¥ 86,000 |
| | 10031228 | qPCR HEX probe assay, 500 reactions | ¥ 48,000 |
| | 10031229 | qPCR HEX probe assay, 1,000 reactions | ¥ 68,000 |
| | 10031230 | qPCR HEX probe assay, 2,500 reactions | ¥ 96,000 |
| | 10031234 | qPCR TEX 615 probe assay, 500 reactions | ¥ 48,000 |
| | 10031235 | qPCR TEX 615 probe assay, 1,000 reactions | ¥ 68,000 |
| | 10031236 | qPCR TEX 615 probe assay, 2,500 reactions | ¥ 96,000 |
| | 10031231 | qPCR Cy5 probe assay, 500 reactions | ¥ 48,000 |
| | 10031232 | qPCR Cy5 probe assay, 1,000 reactions | ¥ 68,000 |
| | 10031233 | qPCR Cy5 probe assay, 2,500 reactions | ¥ 96,000 |
| | 10031237 | qPCR Cy5.5 probe assay, 500 reactions | ¥ 48,000 |
| | 10031238 | qPCR Cy5.5 probe assay, 1,000 reactions | ¥ 68,000 |
| 10031239 | qPCR Cy5.5 probe assay, 2,500 reactions | ¥ 96,000 | |
| mRNA qPCR Probe Assays (Expertly designed — Arabidopsis, chicken, cow, dog, pig, rabbit, rhesus monkey, yeast, zebrafish) | 12001961 | qPCR FAM probe assay, 200 reactions | ¥ 24,000 |
| | 10042982 | qPCR FAM probe assay, 500 reactions | ¥ 34,000 |
| | 10042983 | qPCR FAM probe assay, 1,000 reactions | ¥ 48,000 |
| | 10042984 | qPCR FAM probe assay, 2,500 reactions | ¥ 68,000 |
| | 10042985 | qPCR HEX probe assay, 500 reactions | ¥ 40,000 |
| | 10042986 | qPCR HEX probe assay, 1,000 reactions | ¥ 56,000 |
| | 10042987 | qPCR HEX probe assay, 2,500 reactions | ¥ 78,000 |
| | 10042988 | qPCR TEX 615 probe assay, 500 reactions | ¥ 40,000 |
| | 10042989 | qPCR TEX 615 probe assay, 1,000 reactions | ¥ 56,000 |
| | 10042990 | qPCR TEX 615 probe assay, 2,500 reactions | ¥ 78,000 |
| | 10042991 | qPCR Cy5 probe assay, 500 reactions | ¥ 40,000 |
| | 10042992 | qPCR Cy5 probe assay, 1,000 reactions | ¥ 56,000 |
| | 10042993 | qPCR Cy5 probe assay, 2,500 reactions | ¥ 78,000 |
| | 10042994 | qPCR Cy5.5 probe assay, 500 reactions | ¥ 40,000 |
| | 10042995 | qPCR Cy5.5 probe assay, 1,000 reactions | ¥ 56,000 |
| 10042996 | qPCR Cy5.5 probe assay, 2,500 reactions | ¥ 78,000 | |
| Custom qPCR Probe Assays | 10031261 | Custom qPCR FAM assay, 500 reactions | ¥ 22,000 |
| | 10031262 | Custom qPCR FAM assay, 1,000 reactions | ¥ 32,000 |
| | 10031263 | Custom qPCR FAM assay, 2,500 reactions | ¥ 44,000 |
| | 10031264 | Custom qPCR HEX assay, 500 reactions | ¥ 30,000 |
| | 10031265 | Custom qPCR HEX assay, 1,000 reactions | ¥ 38,000 |
| | 10031266 | Custom qPCR HEX assay, 2,500 reactions | ¥ 52,000 |
| | 10031270 | Custom qPCR TEX 615 assay, 500 reactions | ¥ 34,000 |
| | 10031271 | Custom qPCR TEX 615 assay, 1,000 reactions | ¥ 48,000 |
| | 10031272 | Custom qPCR TEX 615 assay, 2,500 reactions | ¥ 68,000 |
| | 10031267 | Custom qPCR Cy5 assay, 500 reactions | ¥ 34,000 |
| | 10031268 | Custom qPCR Cy5 assay, 1,000 reactions | ¥ 48,000 |
| | 10031269 | Custom qPCR Cy5 assay, 2,500 reactions | ¥ 68,000 |
| | 10031273 | Custom qPCR Cy5.5 assay, 500 reactions | ¥ 34,000 |
| | 10031274 | Custom qPCR Cy5.5 assay, 1,000 reactions | ¥ 48,000 |
| | 10031275 | Custom qPCR Cy5.5 assay, 2,500 reactions | ¥ 68,000 |

Continued

PrimePCR™ Pricing

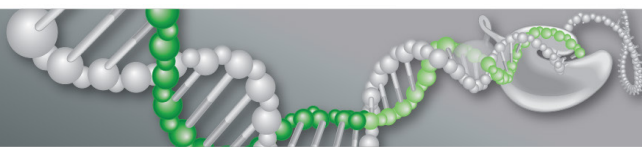


アッセイおよびコントロール

プローブ アッセイ価格

| Product | Catalog # | Description | List Price (JPY) |
|---|-----------|--|------------------|
| lncRNA qPCR Probe Assays (Expertly designed — human) | 12004181 | lncRNA qPCR FAM probe assay, 200 reactions | ¥24,000 |
| | 12004182 | lncRNA qPCR FAM probe assay, 500 reactions | ¥44,000 |
| | 12004183 | lncRNA qPCR FAM probe assay, 1,000 reactions | ¥62,000 |
| | 12004184 | lncRNA qPCR FAM probe assay, 2,500 reactions | ¥86,000 |
| | 12004185 | lncRNA qPCR HEX probe assay, 500 reactions | ¥48,000 |
| | 12004186 | lncRNA qPCR HEX probe assay, 1,000 reactions | ¥67,000 |
| | 12004187 | lncRNA qPCR HEX probe assay, 2,500 reactions | ¥96,000 |
| | 12004188 | lncRNA qPCR TEX 615 probe assay, 500 reactions | ¥48,000 |
| | 12004189 | lncRNA qPCR TEX 615 probe assay, 1,000 reactions | ¥67,000 |
| | 12004190 | lncRNA qPCR TEX 615 probe assay, 2,500 reactions | ¥96,000 |
| | 12004191 | lncRNA qPCR Cy5 probe assay, 500 reactions | ¥48,000 |
| | 12004192 | lncRNA qPCR Cy5 probe assay, 1,000 reactions | ¥67,000 |
| | 12004193 | lncRNA qPCR Cy5 probe assay, 2,500 reactions | ¥96,000 |
| | 12004194 | lncRNA qPCR Cy5.5 probe assay, 500 reactions | ¥48,000 |
| | 12004195 | lncRNA qPCR Cy5.5 probe assay, 1,000 reactions | ¥67,000 |
| | 12004196 | lncRNA qPCR Cy5.5 probe assay, 2,500 reactions | ¥96,000 |

PrimePCR™ Pricing



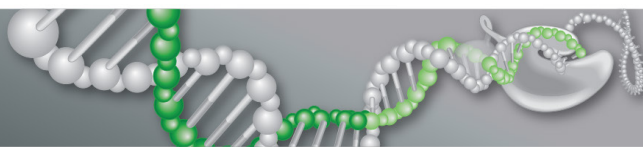
テンプレートおよびコントロールアッセイ

テンプレートおよびコントロールアッセイ価格

| Product | Catalog # | Description | List Price (JPY) |
|---|-----------|---|------------------|
| Positive PCR and Reverse Transcription Control Assays — All Species | 10025591 | PCR control assay, 200 reactions | ¥ 16,000 |
| | 10025695 | RT control assay, 200 reactions | ¥ 16,000 |
| | 10031288 | PCR control FAM assay, 200 reactions | ¥ 20,000 |
| | 10031289 | PCR control HEX assay, 200 reactions | ¥ 20,000 |
| | 10031286 | RT control FAM assay, 200 reactions | ¥ 20,000 |
| Template and Control Assays — Human | 10025716 | Synthetic DNA control template, human | ¥ 16,000 |
| | 10031285 | Synthetic DNA control template probe, human | ¥ 20,000 |
| | 10025694 | RNA quality control assay, human | ¥ 16,000 |
| | 10025352 | gDNA control assay, human | ¥ 16,000 |
| | 10031287 | RNA quality control probe assay, human | ¥ 20,000 |
| | 10031290 | gDNA control probe assay, human | ¥ 20,000 |
| Template and Control Assays — Mouse | 10029101 | Synthetic DNA control template, mouse | ¥ 16,000 |
| | 10031285 | Synthetic DNA control template probe, mouse | ¥ 20,000 |
| | 10029103 | RNA quality control assay, mouse | ¥ 16,000 |
| | 10029105 | gDNA control assay, mouse | ¥ 16,000 |
| | 10032858 | RNA quality control probe assay, mouse | ¥ 20,000 |
| | 10032859 | gDNA control probe assay, mouse | ¥ 20,000 |
| Template and Control Assays — Rat | 10044155 | Synthetic DNA control template, rat | ¥ 16,000 |
| | 10044156 | Synthetic DNA control template probe, rat | ¥ 20,000 |
| | 10044157 | RNA quality control assay, rat | ¥ 16,000 |
| | 10044158 | gDNA control assay, rat | ¥ 16,000 |
| | 10044159 | RNA quality control probe assay, rat | ¥ 20,000 |
| | 10044160 | gDNA control probe assay, rat | ¥ 20,000 |
| Template and Control Assays — Zebrafish | 10047261 | Synthetic DNA control template, zebrafish | ¥ 16,000 |
| | 10047262 | Synthetic DNA control template probe, zebrafish | ¥ 20,000 |
| | 10047263 | RNA quality control assay, zebrafish | ¥ 16,000 |
| | 10047264 | gDNA control assay, zebrafish | ¥ 16,000 |
| | 10047265 | RNA quality control probe assay, zebrafish | ¥ 20,000 |
| | 10047266 | gDNA control probe assay, zebrafish | ¥ 20,000 |
| Template and Control Assays — Yeast | 10047267 | Synthetic DNA control template, yeast | ¥ 16,000 |
| | 10047268 | Synthetic DNA control template probe, yeast | ¥ 20,000 |
| | 10047269 | RNA quality control assay, yeast | ¥ 16,000 |
| | 10047270 | gDNA control assay, yeast | ¥ 16,000 |
| | 10047271 | RNA quality control probe assay, yeast | ¥ 20,000 |
| | 10047272 | gDNA control probe assay, yeast | ¥ 20,000 |
| Template and Control Assays — Rhesus Monkey | 10047273 | Synthetic DNA control template, rhesus monkey | ¥ 16,000 |
| | 10047274 | Synthetic DNA control template probe, rhesus monkey | ¥ 20,000 |
| | 10047275 | RNA quality control assay, rhesus monkey | ¥ 16,000 |
| | 10047276 | gDNA control assay, rhesus monkey | ¥ 16,000 |
| | 10047277 | RNA quality control probe assay, rhesus monkey | ¥ 20,000 |
| | 10047278 | gDNA control probe assay, rhesus monkey | ¥ 20,000 |

Continued

PrimePCR™ Pricing

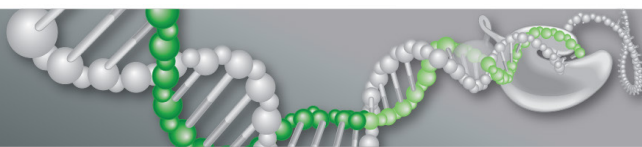


テンプレートおよびコントロールアッセイ

テンプレートおよびコントロールアッセイ価格

| Product | Catalog # | Description | List Price (JPY) |
|--|-----------|--|------------------|
| Template and Control Assays — Pig | 10047279 | Synthetic DNA control template, pig | ¥ 16,000 |
| | 10047280 | Synthetic DNA control template probe, pig | ¥ 20,000 |
| | 10047281 | RNA quality control assay, pig | ¥ 16,000 |
| | 10047282 | gDNA control assay, pig | ¥ 16,000 |
| | 10047283 | RNA quality control probe assay, pig | ¥ 20,000 |
| | 10047284 | gDNA control probe assay, pig | ¥ 20,000 |
| Template and Control Assays — Cow | 10047285 | Synthetic DNA control template, cow | ¥ 16,000 |
| | 10047286 | Synthetic DNA control template probe, cow | ¥ 20,000 |
| | 10047287 | RNA quality control assay, cow | ¥ 16,000 |
| | 10047288 | gDNA control assay, cow | ¥ 16,000 |
| | 10047289 | RNA quality control probe assay, cow | ¥ 20,000 |
| | 10047290 | gDNA control probe assay, cow | ¥ 20,000 |
| Template and Control Assays — Rabbit | 10047315 | Synthetic DNA control template, rabbit | ¥ 16,000 |
| | 10047316 | Synthetic DNA control template probe, rabbit | ¥ 20,000 |
| | 10047317 | RNA quality control assay, rabbit | ¥ 16,000 |
| | 10047318 | gDNA control assay, rabbit | ¥ 16,000 |
| | 10047319 | RNA quality control probe assay, rabbit | ¥ 20,000 |
| | 10047320 | gDNA control probe assay, rabbit | ¥ 20,000 |
| Template and Control Assays — Dog | 10047291 | Synthetic DNA control template, dog | ¥ 16,000 |
| | 10047292 | Synthetic DNA control template probe, dog | ¥ 20,000 |
| | 10047293 | RNA quality control assay, dog | ¥ 16,000 |
| | 10047294 | gDNA control assay, dog | ¥ 16,000 |
| | 10047295 | RNA quality control probe assay, dog | ¥ 20,000 |
| | 10047296 | gDNA control probe assay, dog | ¥ 20,000 |
| Template and Control Assays — Chicken | 10047297 | Synthetic DNA control template, chicken | ¥ 16,000 |
| | 10047298 | Synthetic DNA control template probe, chicken | ¥ 20,000 |
| | 10047299 | RNA quality control assay, chicken | ¥ 16,000 |
| | 10047300 | gDNA control assay, chicken | ¥ 16,000 |
| | 10047301 | RNA quality control probe assay, chicken | ¥ 20,000 |
| | 10047302 | gDNA control probe assay, chicken | ¥ 20,000 |
| Template and Control Assays — <i>Arabidopsis</i> | 10047303 | Synthetic DNA control template, <i>Arabidopsis</i> | ¥ 16,000 |
| | 10047304 | Synthetic DNA control template probe, <i>Arabidopsis</i> | ¥ 20,000 |
| | 10047305 | RNA quality control assay, <i>Arabidopsis</i> | ¥ 16,000 |
| | 10047306 | gDNA control assay, <i>Arabidopsis</i> | ¥ 16,000 |
| | 10047307 | RNA quality control probe assay, <i>Arabidopsis</i> | ¥ 20,000 |
| | 10047308 | gDNA control probe assay, <i>Arabidopsis</i> | ¥ 20,000 |

PrimePCR™ Pricing

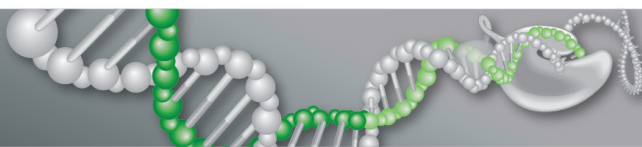


パスウェイおよびコレクションアレイ

アレイ価格

| Product | Catalog # | Description | 最小オーダー 数量 | List Price (JPY) |
|---|---------------------|------------------------|--------------|---------------------|
| 96-Well mRNA qPCR Arrays (SYBR® Green assay) | アレイの種類に よって異なります | 96 well アレイ | | |
| | | 16 assays x 6 samples | 1 | ¥ 26,000 |
| | | 24 assays x 4 samples | 1 | ¥ 26,000 |
| | | 32 assays x 3 samples | 1 | ¥ 30,000 |
| | | 48 assays x 2 samples | 1 | ¥ 32,000 |
| | | 96 assays x 1 sample | 1 | ¥ 34,000 |
| 384-Well mRNA qPCR Arrays (SYBR® Green assay) | アレイの種類に よって異なります | 384 wellアレイ | | |
| | | 16 assays x 24 samples | 1 | ¥ 38,000 |
| | | 24 assays x 16 samples | 1 | ¥ 40,000 |
| | | 32 assays x 12 samples | 1 | ¥ 42,000 |
| | | 48 assays x 8 samples | 1 | ¥ 42,000 |
| | | 96 assays x 4 samples | 1 | ¥ 44,000 |
| | | 128 assays x 3 samples | 1 | ¥ 48,000 |
| | | 192 assays x 2 samples | 1 | ¥ 50,000 |
| | | 384 assays x 1 sample | 1 | ¥ 54,000 |
| 96-Well lncRNA qPCR Arrays (SYBR® Green assay) | アレイの種類に よって異なります | 96 assays x 1 sample | 1 | ¥33,000 |
| 384-Well lncRNA qPCR Arrays (SYBR® Green assay) | アレイの種類に よって異なります | 96 assays x 4 samples | 1 | ¥44,000 |

PrimePCR™ Pricing



カスタムプレート

カスタムプレート価格

| Product | Catalog # | Description | 最小オーダー 数量 | List Price (JPY) |
|---|-----------|----------------------------|--------------|---------------------|
| Custom mRNA 96-Well Plates (SYBR® Green assay) | 10025216 | 96-Well プレート 1-12 genes | 3 | ¥ 24,000 |
| | 10025217 | 13-24 genes | 3 | ¥ 30,000 |
| | 10025218 | 25-36 genes | 3 | ¥ 34,000 |
| | 10025219 | 37-96 genes | 3 | ¥ 40,000 |
| Custom mRNA 384-Well Plates (SYBR® Green assay) | 10025210 | 384-Wellプレート 1-12 genes | 6 | ¥ 34,000 |
| | 10025211 | 13-24 genes | 6 | ¥ 40,000 |
| | 10025213 | 25-36 genes | 6 | ¥ 46,000 |
| | 10025214 | 37-96 genes | 6 | ¥ 52,000 |
| | 10025215 | 97-192 genes | 12 | ¥ 56,000 |
| | 10025212 | 193-384 genes | 12 | ¥ 68,000 |
| Custom lncRNA 96-Well Plates (SYBR® Green assay) | 12004201 | 96-Wellプレート 1-12 genes | 3 | ¥23,000 |
| | 12004202 | 13-24 genes | 3 | ¥29,000 |
| | 12004203 | 25-36 genes | 3 | ¥34,000 |
| | 12004204 | 37-96 genes | 3 | ¥40,000 |
| Custom lncRNA 384-Well Plates (SYBR® Green assay) | 12004205 | 384-Wellプレート 1-12 genes | 6 | ¥34,000 |
| | 12004206 | 13-24 genes | 6 | ¥40,000 |
| | 12004207 | 25-36 genes | 6 | ¥45,000 |
| | 12004208 | 37-96 genes | 6 | ¥51,000 |
| | 12004209 | 97-192 genes | 12 | ¥56,000 |
| | 12004210 | 193-384 genes | 12 | ¥68,000 |

製品の検索、およびご注文の際に必要なオーダーフォームの発行はPrimePCRサイトにて行っていただけます。下記URLよりPrimePCRサイトをご訪問ください。

bio-rad.com/PrimePCR

BIO-RAD バイオ・ラッド ラボラトリーズ 株式会社

ライフサイエンス www.bio-rad.com

本社 〒140-0002 東京都品川区東品川2-2-24 TEL:03-6361-7000 FAX:03-5463-8480

大阪営業所 〒532-0025 大阪市淀川区新北野1-14-11 TEL:06-6308-6568 FAX:06-6308-3064

*学術的お問い合わせは TEL:03-6404-0331 FAX:03-6404-0334

※価格、仕様などは予告無く変更することがありますので、ご了承ください。
 ※価格は2020年11月現在のものです、メーカー希望小売価格（税別）です。
 ※本カタログに記載されている会社名、商品名は各社の商標または登録商標です。

Z12516L 2011a